


RESEARCH ARTICLE

Open Access



Edaravone activates the GDNF/RET neurotrophic signaling pathway and protects mRNA-induced motor neurons from iPSC cells

Qian Li^{1,2}, Yi Feng², Yingchao Xue², Xiping Zhan³, Yi Fu¹, Gege Gui⁴, Weiqiang Zhou⁵, Jean-Philippe Richard⁶, Arens Taga⁶, Pan Li⁵, Xiaobo Mao⁶, Nicholas J. Maragakis⁶ and Mingyao Ying^{2,6*} 

Abstract

Background: Spinal cord motor neurons (MNs) from human iPSCs (iPSCs) have wide applications in disease modeling and therapeutic development for amyotrophic lateral sclerosis (ALS) and other MN-associated neurodegenerative diseases. We need highly efficient MN differentiation strategies for generating iPSC-derived disease models that closely recapitulate the genetic and phenotypic complexity of ALS. An important application of these models is to understand molecular mechanisms of action of FDA-approved ALS drugs that only show modest clinical efficacy. Novel mechanistic insights will help us design optimal therapeutic strategies together with predictive biomarkers to achieve better efficacy.

Methods: We induce efficient MN differentiation from iPSCs in 4 days using synthetic mRNAs coding two transcription factors (Ngn2 and Olig2) with phosphosite modification. These MNs after extensive characterization were applied in electrophysiological and neurotoxicity assays as well as transcriptomic analysis, to study the neuroprotective effect and molecular mechanisms of edaravone, an FDA-approved drug for ALS, for improving its clinical efficacy.

* Correspondence: ying@kennedykrieger.org

²Hugo W. Moser Research Institute at Kennedy Krieger, 707 North Broadway, Baltimore, MD 21205, USA

⁶Department of Neurology, Johns Hopkins University School of Medicine, Baltimore, MD 21205, USA

Full list of author information is available at the end of the article



© The Author(s). 2022 **Open Access** This article is licensed under a Creative Commons Attribution 4.0 International License, which permits use, sharing, adaptation, distribution and reproduction in any medium or format, as long as you give appropriate credit to the original author(s) and the source, provide a link to the Creative Commons licence, and indicate if changes were made. The images or other third party material in this article are included in the article's Creative Commons licence, unless indicated otherwise in a credit line to the material. If material is not included in the article's Creative Commons licence and your intended use is not permitted by statutory regulation or exceeds the permitted use, you will need to obtain permission directly from the copyright holder. To view a copy of this licence, visit <http://creativecommons.org/licenses/by/4.0/>. The Creative Commons Public Domain Dedication waiver (<http://creativecommons.org/publicdomain/zero/1.0/>) applies to the data made available in this article, unless otherwise stated in a credit line to the data.

Results: We generate highly pure and functional mRNA-induced MNs (miMNs) from control and ALS iPSCs, as well as embryonic stem cells. Edaravone alleviates H₂O₂-induced neurotoxicity and electrophysiological dysfunction in miMNs, demonstrating its neuroprotective effect that was also found in the glutamate-induced miMN neurotoxicity model. Guided by the transcriptomic analysis, we show a previously unrecognized effect of edaravone to induce the GDNF receptor RET and the GDNF/RET neurotrophic signaling in vitro and in vivo, suggesting a clinically translatable strategy to activate this key neuroprotective signaling. Notably, edaravone can replace required neurotrophic factors (BDNF and GDNF) to support long-term miMN survival and maturation, further supporting the neurotrophic function of edaravone-activated signaling. Furthermore, we show that edaravone and GDNF combined treatment more effectively protects miMNs from H₂O₂-induced neurotoxicity than single treatment, suggesting a potential combination strategy for ALS treatment.

Conclusions: This study provides methodology to facilitate iPSC differentiation and disease modeling. Our discoveries will facilitate the development of optimal edaravone-based therapies for ALS and potentially other neurodegenerative diseases.

Keywords: Spinal cord motor neuron, Transcription factor, Amyotrophic lateral sclerosis, Synthetic mRNA, MCI-186, Neurotrophic factor

Background

Human pluripotent stem cells (PSCs), including embryonic stem cells (ESCs) and induced pluripotent stem cells (iPSCs), have unique characteristics, such as long-term self-renewal and multi-lineage differentiation capability. Patient-derived iPSCs are widely used in regenerative medicine to provide disease-relevant functional cells for mechanistic studies, drug discovery and cell replacement therapy [1–3]. All iPSC-based applications rely on robust differentiation strategies for the manufacture of lineage-specific and functional progenies.

Human motor neurons (MNs) derived from iPSCs provide a unique and efficient platform for modeling various MN disorders and developing effective therapies. Amyotrophic lateral sclerosis (ALS) is a rapidly progressive neurodegenerative disease that is characterized primarily by MN degeneration in the brain and spinal cord and has a median survival of 20 to 48 months [4]. Therapeutic development for ALS is extremely challenging [5]. The Food and Drug Administration (FDA) has only approved two drugs with modest efficacy. The first FDA-approved ALS drug Riluzole is a glutamatergic neurotransmission inhibitor, and only improves patient survival by 2–3 months without showing benefit on motor function [6–8]. The recently approved edaravone (also known as Radicava or MCI-186) was shown to slow early-stage disease progression in a subset of ALS patients enrolled in a phase III study [9]. However, controversy remains over the clinical efficacy of edaravone in ALS patients [10–12]. It is also unclear how edaravone might be effective in ALS, although it is predicted to act through reducing oxidative stress based on its proposed function as a free radical scavenger [13, 14]. To improve clinical efficacy of current ALS drugs, it is critical to understand their molecular mechanisms of action in

disease models that closely recapitulate the genetic and phenotypic complexity of ALS, and further rationally design single or possibly combination therapies together with predictive biomarkers to achieve better clinical efficacy.

ALS is a highly heterogeneous disease with regard to phenotype and progression, a large number of disease-associated genetic variants, and multiple cellular pathways affected [5]. A bank of iPSC-derived MNs from familial and sporadic ALS patients provides an efficient modeling system mimicking the complexity of ALS and suitable for high-throughput drug screening. Building this MN bank needs a robust iPSC differentiation strategy. Traditional MN differentiation methods mainly use small molecule compounds to induce neural conversion and MN lineage specification in iPSCs. These multi-step protocols take about 10–14 days to generate MN precursors with variable purity (commonly 50–70%) [15, 16]. Since the principle of all MN differentiation strategies is based on generating Olig2⁺ MN progenitors [15], it may be feasible to drive robust MN conversion through ectopic expression of Olig2 and other transcription factor (TF) drivers of MN development (e.g. Ngn2) [17, 18]. Ectopic TF expression can be achieved by synthetic mRNA delivery, an efficient, non-viral and non-integrating strategy that has been used by us to differentiate iPSCs to dopaminergic neurons [19].

Here, we established a rapid MN differentiation method using synthetic mRNAs coding two TFs. These MNs were applied to high-throughput phenotypic analysis and transcriptomic profiling to study the ALS drug edaravone. We revealed a novel effect of edaravone in activating the neurotrophic factor signaling pathway in vitro and in vivo, and found an effective combination strategy to protect MNs.

Materials and methods

Cell culture

The control iPSC line N1 and N3 (referred to as iPSC1 and iPSC3) was derived from human skin fibroblasts as previously characterized and used by us [19, 20]. Two ALS iPSC lines (referred to as iPSC2 and iPSC4) was derived from skin fibroblasts from an ALS patient with the *SOD1*^{A4V} mutation. Cell reprogramming was performed using the Sendai virus system with SOX2/OCT4/KLF4/MYC (CytoTune-iPS Reprogramming Kit; Thermo Fisher Scientific, Rockville, MD, <http://www.thermofisher.com>). iPSC pluripotency is characterized by immunocytochemistry for pluripotent markers (NANOG, OCT4, TRA-1-60, and SSEA-3) and embryoid body formation assay. The human ESC line H1 was obtained from WiCell Research Resources (Madison, WI, <http://www.wicell.org>). iPSCs and ESCs were maintained in mTESR1 medium (Stem Cell Technologies, Vancouver, BC, Canada) at 5% CO₂/95% air condition at 37 °C and were passaged using ReLeSR™ (Stem Cell Technologies). Karyotype analysis of G-banded metaphase chromosomes has been performed to confirm chromosomal integrity. Human neural stem cells were established from a fetal brain, immortalized by *v-myc* and extensively characterized by Dr. Vescovi and his colleagues [21]. Astrocyte differentiation uses DMEM/F12 medium (Thermo Fisher Scientific) with 1% fetal bovine serum (FBS) for 21 days following the publication [21].

All chemicals were from Sigma Aldrich unless otherwise mentioned.

mRNA synthesis and transfection

Coding sequences of human Ngn2 and Olig2 were cloned into a vector containing the T7 promoter and poly(A) tail for in-vitro transcription as reported by us [19]. mRNA transfection used the Lipofectamine™ Stem Transfection Reagent (Thermo Fisher Scientific). For each well of the 12-well plate, we used 0.25 µg mRNAs (N-SA:O-SA = 1:1) with 1.5 µl lipid. All procedures involving recombinant DNA follow the National Institutes of Health guidelines.

mRNA-induced MN differentiation

iPSCs were plated at a density of 3×10^5 cells per well in a 12-well plate coated with growth-factor-reduced Matrigel (Corning). iPSCs were transfected daily with N-SA/O-SA mRNAs for 3 days. Culture medium with SHH (100 ng/ml) and DAPT (10 µM) were changed daily and shifted from mTeSR1 to N2 (Thermo Fisher Scientific) in 3 days. Cells were dissociated by Accutase and replated to poly-D-Lysine/Laminin-coated surface at the density of 1×10^5 cells/cm². Neuron maturation medium contain neurobasal medium with the B27 supplement, BDNF (10 ng/ml), GDNF (10 ng/ml), cAMP (0.1 mM),

ascorbic acid (0.2 mM), DAPT (10 µM). Medium was changed after 48 h followed by half change every 3–4 days. Cryopreservation medium containing 40% neurobasal medium, 50% FBS and 10% DMSO.

Western blotting

We followed our previous publication [19] to perform total protein extraction using RIPA buffer (Sigma-Aldrich), SDS-polyacrylamide gel electrophoresis and western blotting. Protein levels were quantified with the Odyssey IR Imaging System (LI-COR Biosciences). All primary antibodies are listed in Table S6. The original blot images were included in the additional file “western_blot.pdf”.

Immunofluorescence staining and quantification

Cells were fixed in 4% paraformaldehyde in PBS (pH 7.4) and subjected to immunostaining as published by us [19], using primary antibodies listed in Table S6. The percentage of marker-positive cells over DAPI+ nuclei was determined in samples from at least three samples that were independently differentiated for staining. The high-content analysis software (HCA-Vision V2.2.0. CSIRO) was used for nucleus detection and cell body segmentation. Threshold for each marker was set based on signal intensity in the IgG isotype control.

GDNF ELISA

ELISA used the kit from Thermo Fisher Scientific (human GDNF) and Rockland (mouse GDNF).

Electrophysiological recordings

Voltage-clamp recording was performed at 35 °C in a chamber perfused with regular artificial cerebrospinal fluid (124 mM NaCl, 2.5 mM KCl, 1.3 mM MgCl₂, 2.5 mM CaCl₂, 1 mM NaH₂PO₄, 26.2 mM NaHCO₃, 20 mM glucose, pH 7.4, equilibrated with 95% O₂ and 5% CO₂, ~310 mosm), which flowed at 3 ml/minute. Patch electrodes were pulled from borosilicate glass and had resistances of 2.0–4.0 MΩ when filled with an intracellular solution (135 mM KMeSO₄, 5 mM KCl, 5 mM HEPES, 0.25 mM EGTA-free acid, 2 mM Mg-ATP, 0.5 mM GTP, 10 mM phosphocreatine-tris, pH 7.3, ~290 mosm).

Neurons were identified using a 10X objective mounted on an upright microscope with transmitted light, and their neuronal somata were then visualized through a 40X water immersion objective using IR differential interference contrast optics (DIC). The cell somatic recordings were made using an Axopatch 700B amplifier in combination with pClamp 11 software (Molecular Devices). Neurons were initially voltage-clamped at –70 mV, and R_{series} and R_{input} were monitored using a 2.5-mV 100-ms depolarizing voltage step in each recording sweep. The current traces were filtered at 5 kHz,

digitized at 10 kHz using a Digidata 1550b interface, and stored for off-line analysis. Next, recording was switched to current clamp. The resting membrane potential and the action potential were monitored for more than 5 min before drug applications. To induce action potentials, the neurons were commanded by multiple steps of hyperpolarization currents. Tetraethylammonium (TEA-Cl, 1 mM) and TTX (0.5 μ M) from Sigma-Aldrich were added to the artificial cerebrospinal fluid, to block K⁺ or Na⁺ channels, respectively. Electrophysiological recording data were first visualized with Clampfit 11 and exported to MATLAB (Mathworks, Natick, MA, <http://www.mathworks.com>) for analysis. The recording traces were visualized using Igor Pro 6.0 (WaveMetrics, Portland, OR, <http://www.wavemetrics.com>).

Multi-electrode Array (MEA)

After 4 days of differentiation, neurons were plated on poly-D-lysine/laminin coated CytoView MEA 24-well plates (Axion BioSystems, www.axionbiosystems.com) at a density of 5×10^4 cells/well. Recordings from electrodes were made using a Maestro MEA system (Axion BioSystems). Data were sampled at 12.5 kHz, digitized, and analyzed using the Axion Integrated Studio software (Axion BioSystems) with a 200 Hz high pass and 4 kHz low pass filter and an adaptive spike detection threshold set at 6 times the standard deviation of the background noise for each electrode with 1 s binning. Total action potential counts, mean neuronal firing rates and total burst counts were quantified using the Axion Integrated Studio software.

For H₂O₂ and edaravone treatment, a 10-min MEA recording was acquired as the initial baseline at 0 h. Neurons were treated with edaravone followed by adding H₂O₂ to the edaravone-containing medium. Spontaneous action potential parameters after H₂O₂ treatment were normalized to the initial baseline.

Neurotoxicity assay, Calcein-AM staining, neurite tracing and high-content analysis

Neurons (1×10^3 per well) were plated in plates coated with poly-D-lysine/laminin with 10^3 and 10^4 per well for the 1536- and 96-well plate, respectively. Neurons were pre-treated with edaravone (10 μ M) for 16 h in neurotrophic factor-free neurobasal medium followed by H₂O₂ (25 μ M) or glutamate (200 μ M) treatment for 24 h. Live cell imaging used Calcein-AM dye (1 μ M, Thermo Fisher Scientific). Neurite length quantification used the high-content analysis software (HCA-Vision V2.2.0, CSIRO). Neurite length per field was normalized to the number of Hoechst 33342 stained nuclei.

RNA sequencing and quantitative real-time PCR (qPCR)

Total RNAs were extracted using the RNeasy Mini kit (Qiagen) and subjected to sequencing using the HiSeq 2500 platform (Illumina). Raw reads were aligned to reference human genome build hg19 using HISAT2 [22] with default parameters. For each gene, the number of reads aligned to its exons were counted and summarized into gene level counts by StringTie [23] based on the GENCODE hg19 annotation. Normalization between samples was carried out by the R package edgeR [24, 25], which controls sequencing depth and RNA composition effects. Heatmap was generated according to the count table with scaling across the samples for each gene. The RNA-Seq data sets can be accessed through the Gene Expression Omnibus (GEO) Repository (GSE151997).

qPCR analysis follow our previous publication [19]. Relative expression of each gene was normalized to the 18S rRNA. Primer sequences are listed in Table S6.

Edaravone treatment in mice

C57BL/6J mice (Jackson Laboratory, 8 weeks old, female) received daily intraperitoneal administration of vehicle (saline) or edaravone (15 mg/kg body weight), following the previous publication [26]. Entire spinal cord tissues were harvested using the hydraulic extrusion method [27]. Total proteins were extracted using RIPA buffer containing protease and phosphatase inhibitors.

Availability of data and materials

Further information and requests for resources and reagents will be fulfilled by the corresponding author. All unique/stable reagents generated in this study are available with a completed Materials Transfer Agreement. The RNA-Seq datasets reported here have been deposited to the Gene Expression Omnibus (GEO) Repository (GSE151997).

Study approval

All experiments involving human stem cells have been approved by the Johns Hopkins Medicine Institutional Review Board. The animal protocol was approved by the Johns Hopkins School of Medicine Animal Care and Use Committee.

Statistics

All quantifications were performed by observers blinded to the experimental groups. All results represent at least three replicates with details in each figure legend. All data are represented as Mean \pm SEM. Statistical analysis was performed using Prism software (GraphPad, San Diego, CA, <http://www.graphpad.com>). For comparing two groups, unpaired, two-tailed Student's t test was

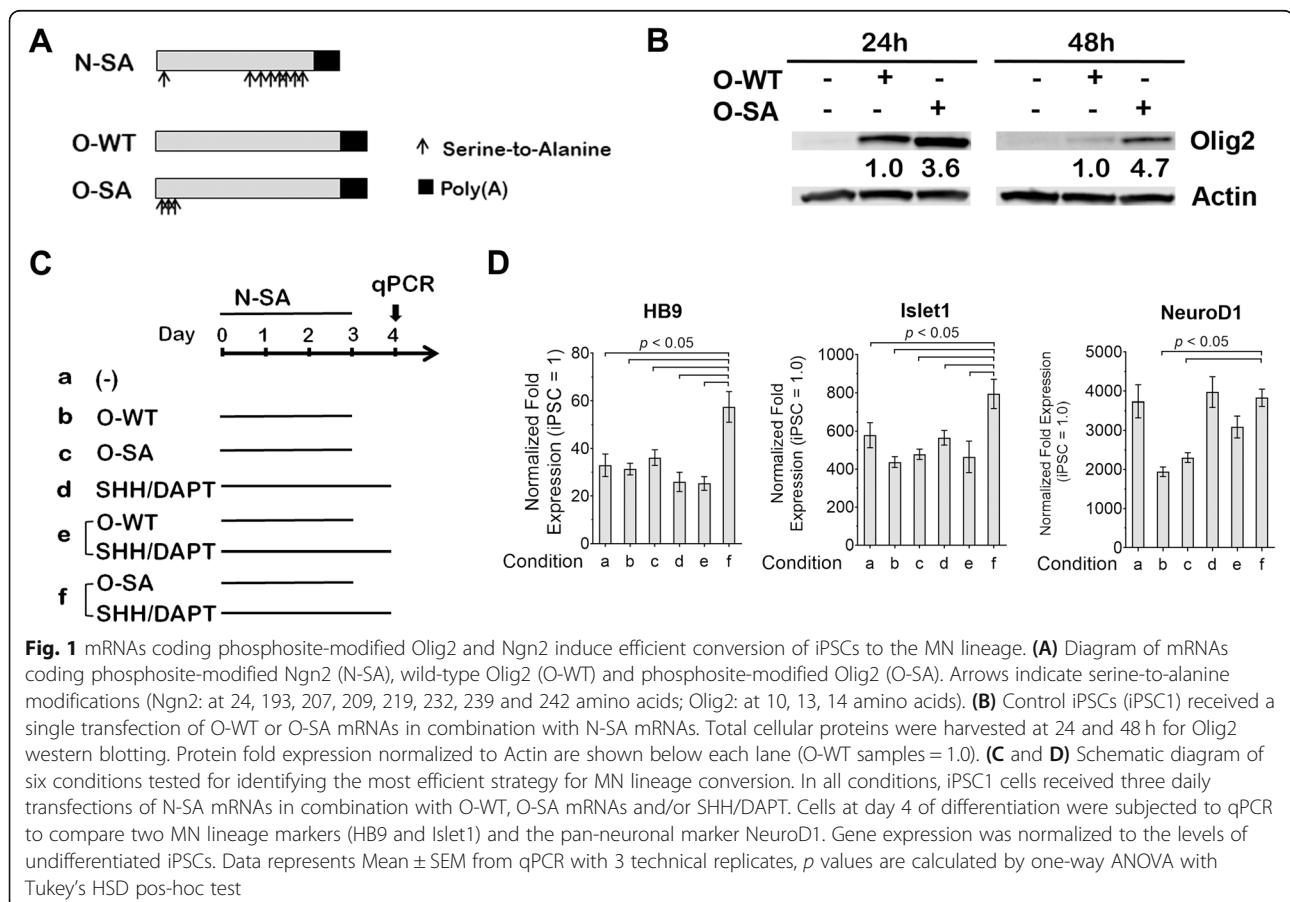
performed (minimal requirement: $p < 0.05$). For more than two groups, one-way ANOVA with Tukey's honestly significant difference (HSD) post-hoc test was used (minimal requirement: $p < 0.05$). qRT-PCR analysis used t test with the Bonferroni correction for multiple comparisons. Other statistical tests were specified in each figure legend.

Results

Synthetic mRNAs coding Olig2 and Ngn2 with phosphosite modification induce efficient MN differentiation from iPSCs

Ngn2 and Olig2 are two TFs co-expressing in motor neuron progenitor cells, and their ectopic expression in combination induces MNs in the chick neural tube [17]. Intrigued by these results, our goal is to develop synthetic mRNAs to ectopically express Ngn2 and Olig2 in iPSCs for efficient generation of mRNA-induced MNs, hereinafter referred to as miMNs. We previously reported that mRNAs coding Ngn2 with eight serine-to-alanine modifications (Fig. 1A, referred to as N-SA) lead to higher protein expression and more efficient neuronal conversion in iPSCs, compared to mRNAs coding wild-type Ngn2 [19]. Thus, we use N-SA mRNAs here for

MN differentiation. To optimize Olig2 mRNAs, we compared mRNAs coding wild-type Olig2 and a modified form with three serine-to-alanine mutations at S10, S13 and S14 sites (Fig. 1A, referred to as O-WT and O-SA, respectively). We co-transfected iPSCs using N-SA mRNAs in combination with two forms of Olig2 mRNAs to determine Olig2 protein levels in this context with Ngn2 expression. O-SA mRNAs produced 2.6-fold and 3.7-fold more proteins than O-WT mRNAs, at 24 h and 48 h after transfection, respectively (Fig. 1B), suggesting that O-SA mRNAs may more efficiently induce MN differentiation. We further tested six differentiation schemes (Fig. 1C) that combine N-SA with O-WT or O-SA mRNAs, also including SHH, a known morphogen for MN lineage specification [28], and DAPT, a Notch signaling inhibitor widely used to promote neuronal conversion [29]. Three daily co-transfections of N-SA/O-SA mRNAs plus SHH/DAPT most efficiently induced MN lineage conversion, based on significantly higher levels of HB9 and Islet1, two well-defined MN lineage markers (Fig. 1D). This combination is also among the top-three schemes showing equal efficiency in inducing the neuronal marker NeuroD1 (Fig. 1D). In contrast, other schemes (e.g. N-SA/O-WT/SHH/DAPT) did not

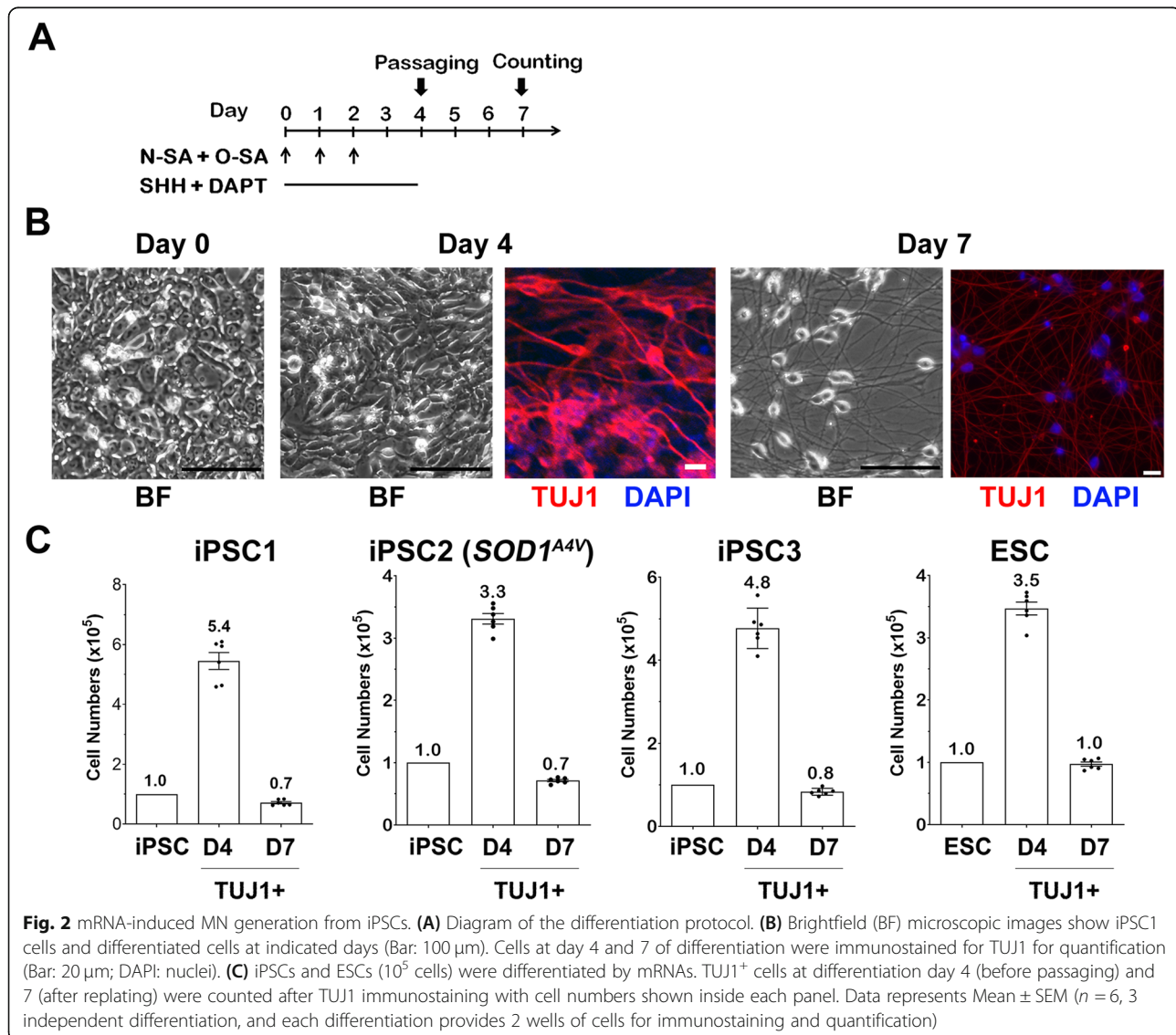


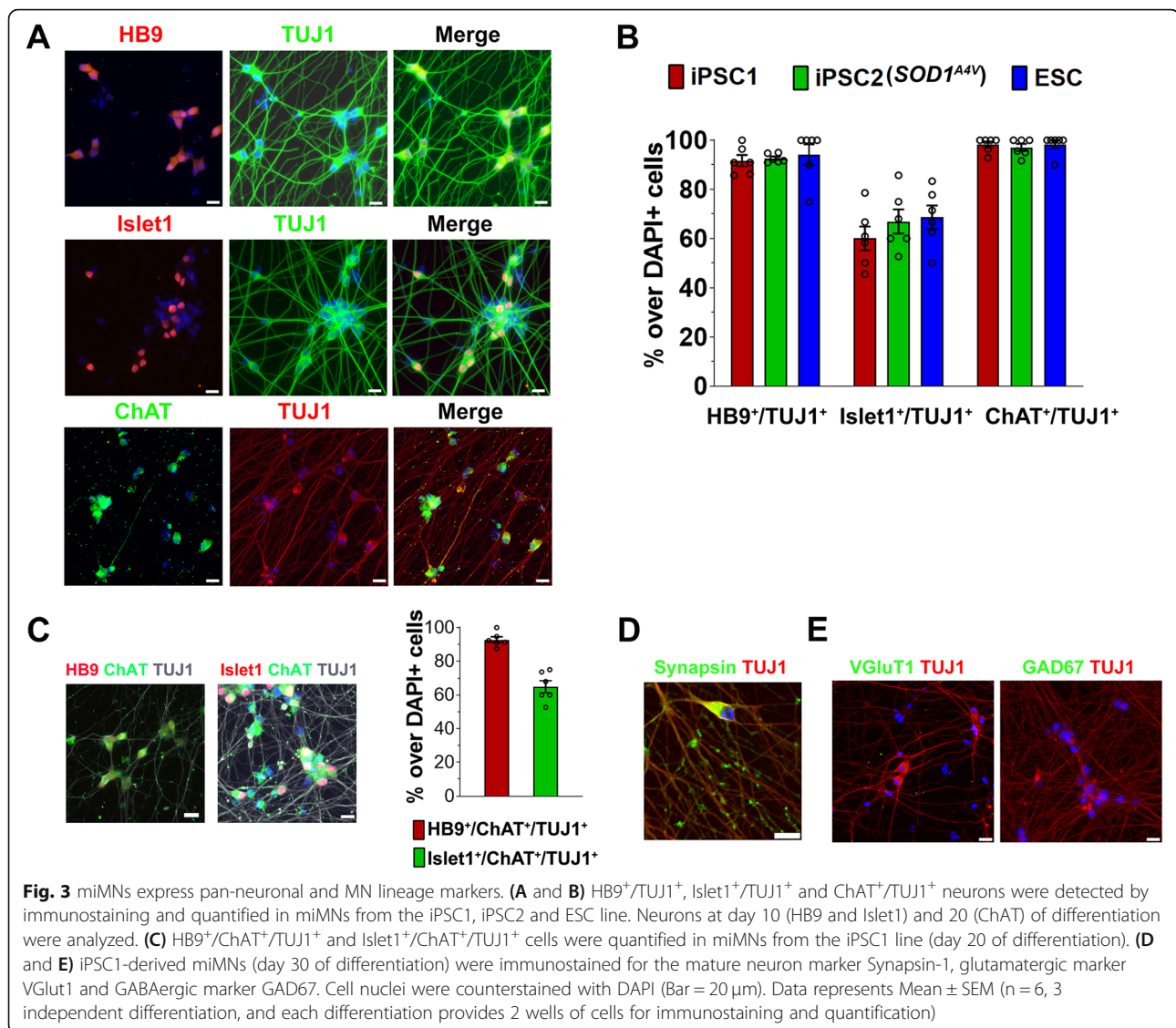
show similar efficiency in simultaneously inducing neuronal and MN lineage markers.

Guided by these results, we established a rapid 4-day MN differentiation protocol, including three daily co-transfections of N-SA/O-SA mRNAs plus SHH/DAPT treatment (Fig. 2A, also see Materials and Methods). In 4 days, mRNA transfection converted iPSCs to TUJ1⁺ neuronal cells (Fig. 2B, TUJ1⁺: >90%). These cells, after being passaged, can be cryopreserved or matured in vitro to miMNs showing typical neuronal morphology (Fig. 2B, TUJ1⁺: >95%) and expressing various MN markers as characterized below. This protocol reproducibly generated miMNs from two control iPSC lines (iPSC1 and iPSC3), an ALS iPSC line with the SOD1^{A4V} mutation (iPSC2), as well as human ESCs (Fig. 2C).

miMNs from control and ALS iPSCs as well as ESCs express known MN markers, including HB9, Islet1 and

the cholinergic neuron marker ChAT (Fig. 3A and B; HB9⁺/TUJ1⁺: >92%; Islet1⁺/TUJ1⁺: >70%; ChAT⁺/TUJ1⁺: >97%). miMNs also co-express MN markers (Fig. 3C; HB9⁺/ChAT⁺/TUJ1⁺: >93%; Islet1⁺/ChAT⁺/TUJ1⁺: >65%). The pluripotent stem cell marker (OCT4) and the oligodendrocyte lineage marker (O4) were not detected in miMNs (Supplemental Fig. 1A). The cholinergic neuron marker ChAT is more abundantly expressed by miMNs at day 20 of differentiation (Fig. 3A and B), compared to those at day 10 of differentiation (Supplemental Fig. 1B). miMNs after in-vitro maturation also express the synaptic vesicle protein and mature neuron marker Synapsin 1 along TUJ1⁺ nerve fibers (Fig. 3D). Ngn2 induction alone has been reported to induce non-MN subtypes in iPSCs and ESCs (e.g. cortical glutamatergic neurons) [30, 31]. Thus, we assessed markers for glutamatergic (VGluT1) and GABAergic





(GAD67) neurons and showed the absence of these markers in miMNs (Fig. 3E), supporting MN conversion driven by N-SA and O-SA in combination. As a control, iPSCs were differentiated by three daily transfections of NSA mRNA alone without using OSA mRNA or SHH/DAPT. NSA-induced TUJ1⁺ neurons express the glutamatergic marker VGLut1 (Supplemental Fig. 1C), which result is consistent with publications showing glutamatergic neuron differentiation driven by Ngn2 alone [30, 31].

Electrophysiological and high-content analysis of miMNs

The functional maturation of miMNs was studied using patch-clamp recording (Fig. 4A). In miMNs from the control iPSC line iPSC1, all of 36 recorded neurons (15–20 days of in-vitro maturation) showed spiking following steps of hyperpolarization current injections (Fig. 4B).

Repetitive multiple action potentials were induced by the hyperpolarization in 19 of 36 neurons (Fig. 4C, left panel). The action potentials were confirmed by blocking with TTX, a selective sodium channel blocker [32] (Fig. 4C). As the generation of action potentials requires a coordinated interplay between sodium and potassium channels, we showed that TEA, a voltage-dependent K⁺ channel blocker [33] attenuated the repetitive action potentials that recovered after compound washing-out (Fig. 4D). In addition, the late interspike intervals (ISIs) were significantly longer than that for the preceding ISIs (Fig. 4E; 24.3 ± 1.3 vs 13.8 ± 1.2 ms, $p < 0.001$), implicating that miMNs exhibit adaptation to extending hyperpolarization, a previously reported physiological property of bona fide MNs [34–36].

Next, we applied the microelectrode array (MEA) system to continuously monitor spontaneous neuronal

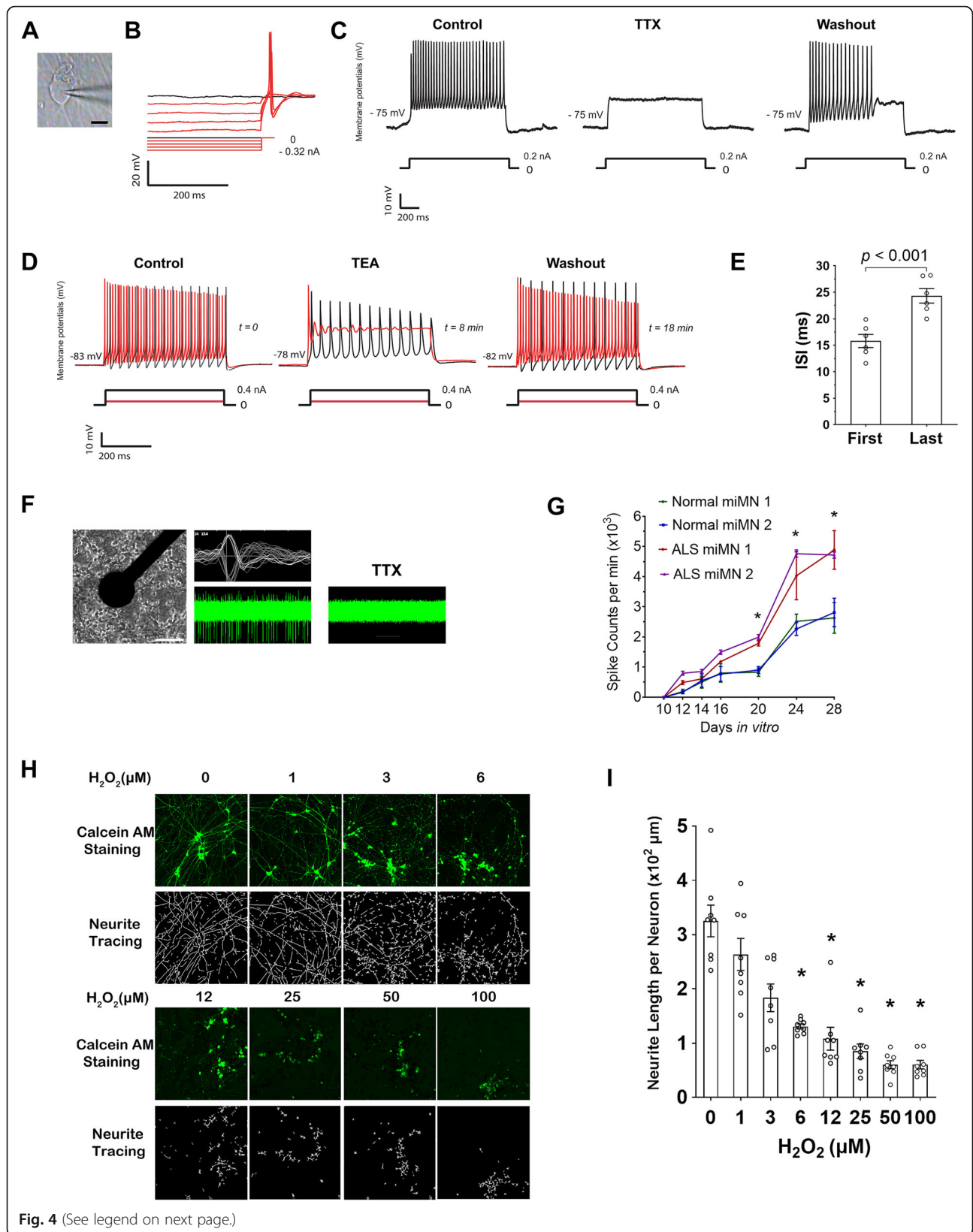


Fig. 4 (See legend on next page.)

(See figure on previous page.)

Fig. 4 Functional characterization and high-content analysis of miMNs. **(A)** Microscopic image of a patched miMN (Bar = 10 μ m). **(B)** Typical rebound potentials after steps of hyperpolarization in miMNs. **(C and D)** miMNs displayed repetitive action potentials after depolarization commands. The action potentials were reversibly blocked by TTX (C) and TEA (D). Traces underneath each panel indicate current steps. **(E)** Interspike intervals (ISI) to depolarization current injections in miMNs are significantly delayed ($n = 6$, $p = 0.001$, t-test). The first ISI indicates the ISI between the first and second spike following depolarization commands, whereas the last ISI indicates the ISI between the last two spikes after depolarization commands being withdrawn. **(F)** A brightfield image shows miMNs attached on the electrode underneath the MEA plate (Bar = 100 μ m). The right panels show spiking waveforms (top) recorded by one electrode and spontaneous spiking activity of miMNs (bottom) with +/- TTX treatment (0.5 μ M). **(G)** Control miMNs from iPSC1 and iPSC3 iPSCs (referred to as control miMN 1 and 2, respectively) and ALS miMNs from iPSC2 and iPSC4 iPSCs (referred to as ALS miMN 1 and 2, respectively), were plated to the MEA plate at day 4 of differentiation. Their spontaneous spiking was recorded at indicated days in vitro. ALS miMNs showed more spiking activity than control miMNs (3 technical replicates for each miMN line, *: $p < 0.05$, linear regression with clustered data, ALS vs control miMNs). **(H and I)** miMNs from iPSC1 at day 4 of differentiation were cultured in the 1536-well plate for 2 days, and subjected to 24 h H_2O_2 treatment followed by Calcein AM staining, neurite tracing (H) and neurite length quantification (I). Neurite length was normalized to nuclei numbers (6 wells for each condition as technical replicates, *: $p < 0.01$, one-way ANOVA with Tukey's HSD pos-hoc test, compared to the untreated control). Data represents Mean \pm SEM.

activities of miMNs in a high-throughput manner. miMNs formed direct contact with electrodes for recording spontaneous spikes that were blocked by the selective sodium channel blocker TTX (Fig. 4F). During in-vitro maturation, control miMNs from two iPSC lines showed more active spontaneous firing and their spiking rate reached the peak after 24 days (Fig. 4G). ALS miMNs from two iPSC lines with the *SOD1*^{A4V} mutation showed more active spontaneous firing than control miMNs (Fig. 4G, $p < 0.05$), consistent with the previous report showing hyperexcitability of MNs from ALS iPSCs [37]. The number of active electrodes did not show difference between control and ALS miMNs (Supplemental Fig. 1D). At the end of the MEA analysis (30 days in vitro), neuron numbers in MEA wells showed no difference between control and ALS miMNs ($0.8\text{--}1.3 \times 10^4$ per well).

miMNs are applicable to high-content analysis in the 1536-well format. miMNs from iPSC1 at day 4 of differentiation were plated using a microplate dispenser. After 48 h, miMNs stained with Calcein-AM showed robust neurite outgrowth (Fig. 4H). These miMNs were treated with hydrogen peroxide (H_2O_2) that has been used to model oxidative stress and neurotoxicity in ALS and other neurodegenerative diseases [38]. H_2O_2 dose-dependently induced morphological signs of neurotoxicity, including neurite fragmentation and cell body condensation (Fig. 4H and I, $IC_{50} = 5.4 \mu$ M, $p < 0.01$, compared to the untreated control).

Edaravone protects miMNs from H_2O_2 -induced neurotoxicity

We applied miMNs to explore the cellular and molecular effects of edaravone, an FDA-approved drug for ALS. In H_2O_2 -induced neurotoxicity assay, edaravone significantly alleviated neurite damage in miMNs (Fig. 5A and B). Edaravone-treated miMNs showed only 26% reduction of neurite length after H_2O_2 (25 μ M) treatment, compared to the edaravone-untreated control ($p < 0.05$).

This neurite damage was less severe than the 93% reduction of neurite length shown in edaravone-untreated neurons ($p < 0.01$). Without H_2O_2 treatment, edaravone did not significantly alter neurite length in miMNs. The neuroprotective effect of edaravone was further tested using the MEA system. We used H_2O_2 at a low concentration (3 μ M) that does not significantly damage neurites as shown in Fig. 4H and I). This low-dose H_2O_2 treatment still inhibited spontaneous spiking of miMNs by about 70% (Fig. 5C and D), recapitulating the early impairment of neuronal functions before neurodegeneration. In H_2O_2 -treated miMNs, edaravone effectively restored neuronal spiking to the same level as H_2O_2 -untreated neurons (Fig. 5C and D, $p < 0.05$). In H_2O_2 -untreated miMNs, edaravone did not significantly alter spontaneous spiking, indicating that this drug may function through neuroprotection but not through elevating neuronal activity.

We also studied glutamate-induced neurotoxicity in miMNs and found that edaravone significantly alleviated glutamate-induced neurite damage in miMNs (Fig. 5E and F). Edaravone-treated miMNs showed only a 15% reduction of neurite length after glutamate treatment (200 μ M, 24 h), significantly less than the 57% reduction in edaravone untreated neurons ($p < 0.05$). Taken together, these results support the neuroprotective function of edaravone in miMNs, thus warranting studies to reveal underlying mechanisms.

Transcriptomic profiling of edaravone-induced molecular responses in miMNs

We performed RNA-sequencing to identify differentially-expressed (DE) genes and their associated signaling pathways. The control iPSC line (iPSC1) was independently differentiated into two sets of miMNs as technical replicates. After in-vitro maturation for 20 days, these two sets of miMNs were treated with edaravone (10 μ M) or DMSO as the control for 24 h. This edaravone concentration has been shown to protect

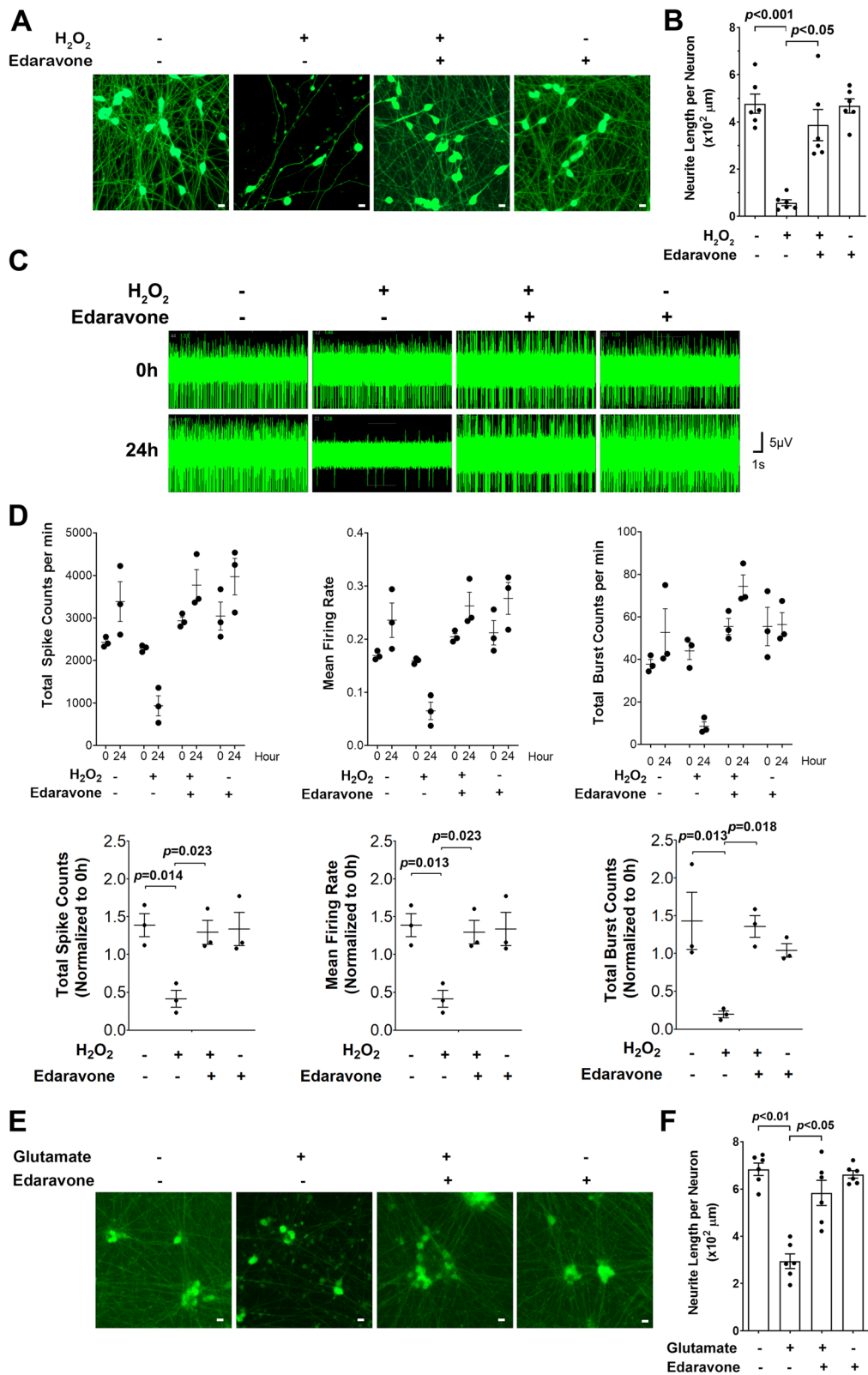


Fig. 5 (See legend on next page.)

(See figure on previous page.)

Fig. 5 Edaravone protects miMNs from H₂O₂-induced neurotoxicity. **(A and B)** iPSC1-derived miMNs (day 7 of differentiation) were pre-treated with edaravone (10 μM) for 16 h in neurotrophic factor-free medium followed by H₂O₂ treatment (25 μM, 24 h). After Calcium AM staining (A, Bar = 10 μm), neurite length was quantified (B, 6 wells for each condition as technical replicates). Edaravone rescues H₂O₂-induced neurite damage. **(C and D)** iPSC1-derived miMNs (day 25 of differentiation) were pre-treated with edaravone (10 μM) in neurotrophic factor-free medium for 16 h followed by H₂O₂ treatment (3 μM, 24 h). Representative traces **(C)** show spontaneous spiking activity that was quantified **(D, top panel)** and compared using values normalized to 0 h **(D, bottom panel)**. Each condition has 3 wells as technical replicates. Edaravone restores spontaneous spiking activity impaired by H₂O₂. **(E and F)** iPSC1-derived miMNs (day 7 of differentiation) were pre-treated with edaravone (10 μM) for 16 h in neurotrophic factor-free medium followed by glutamate treatment (200 μM, 24 h). After Calcium AM staining (E, Bar = 10 μm), neurite length was quantified (F, 6 wells for each condition as technical replicates). Edaravone rescues glutamate-induced neurite damage. Data represents Mean ± SEM with *p* values (one-way ANOVA with Tukey's HSD pos-hoc test) shown inside each panel

miMNs from H₂O₂- and glutamate-induced neurotoxicity (Fig. 5). Over 20 million cDNA reads were generated for each of the two conditions (two technical replicates for each condition) and showed a more than 90% alignment rate to the human genome. Consistency between two replicates was demonstrated by heatmap clustering (Supplemental Fig. 2A). We also performed transcriptomic comparison between miMNs and MNs differentiated by the traditional compound-based method [39]. miMN samples more closely cluster with compound-induced MNs, especially those at 8–9 days of differentiation (Supplemental Fig. 2B), compared to cells at the undifferentiated (Day 0) and early-induction (Day 1–2) stage.

There were 2329 up-regulated (Table S1) and 1916 down-regulated genes (Table S2) altered by edaravone (Fig. 6A, FDR ≤ 0.01, log₂(Fold-Change) Cut-off = ±0.5, see Table S3 for expression data of all detected genes). The Ingenuity Pathway Analysis (IPA) was used for gene functional annotation and pathway enrichment assay. IPA pathways enriched in up- and down-regulated genes were ranked in Table S4 and S5, respectively. Top ten IPA pathways (Fig. 6B) include those associated with neuron functions and ALS pathogenesis, such as the synaptogenesis and CREB signaling enriched in up-regulated genes, and the mitochondrial dysfunction and oxidative phosphorylation signaling enriched in down-regulated genes. Four IPA pathways and associated DE genes were chosen for further studies (Fig. 6C with highlighted genes for validation), including the ALS, superoxide radical degradation, GDNF and neurotrophin/TRK signaling pathways. This selection was based on essential roles of these pathways and their associated genes in MN survival and function, and ALS pathogenesis.

We used qPCR to validate DE genes of interest in both control and ALS miMNs (Fig. 6D and E). Validated up-regulated genes include key components of the neurotrophic factor signaling, such as *GDNF* and its receptor *RET*, another neurotrophic gene *VGF*, as well as two downstream signaling molecules (*GRB2* and *RASD2*). The GDNF co-receptor *GFRA1* is a down-regulated gene in the RNA-seq result, whereas qPCR did not detect its

change in two miMN models. Validated up-regulated genes also include two antioxidant enzymes (*CAT* and *GPX7*). The up-regulation of some ALS signaling-associated genes was validated in control and ALS miMNs, including *GRID2* and *GRIK4* from the glutamate receptor family. *GRIA1* upregulation is validated in control miMNs. Down-regulated genes that were consistently validated in control and ALS miMNs include the pro-apoptotic gene *BAX* and *CYGB* coding a stress-responsive hemoprotein expressed in the brain [40]. *SOD1* showed a small (15%) but significantly down-regulation in control but not ALS miMNs. *SOD2* showed about 40% down-regulation only in ALS miMNs.

Edaravone promotes the GDNF/RET neurotrophic signaling pathway in miMNs and the mouse spinal cord

Consistent with qPCR results, edaravone elevated the levels of RET and VGF proteins from the neurotrophic factor signaling and two antioxidant enzymes (*CAT* and *GPX7*), in control and ALS miMNs (Fig. 7A). Protein levels of the GDNF co-receptor GFRA1 were also elevated by edaravone in both miMN models (Fig. 7A), although GFRA1 mRNAs were not altered by edaravone as previously determined by qPCR (Fig. 6D and E). Next, we focused on the previously unrecognized effect of edaravone on the GDNF/RET neurotrophic signaling. Edaravone treatment in control and ALS miMNs induces the expression of HOXB5, NKX2.1 and PHOX2B (Fig. 7B), three TFs transactivating the RET promoter [41, 42], suggesting the involvement of these TFs in activating RET transcription by edaravone. As a control, we tested other TFs (CTCF, MYC, RAD21 and GABPA), the binding motifs of which are found in the RET promoter. We did not detect their induction by edaravone (data not shown). To determine the functional outcome of GDNF receptor induction by edaravone, we performed GDNF stimulation (30 min) in control and ALS miMNs that have been cultured overnight in neurotrophic factor-free medium with or without edaravone. Edaravone-treated miMNs showed higher levels of total and phosphorylated RET proteins than the untreated control (Fig. 7C). More importantly, edaravone-treated

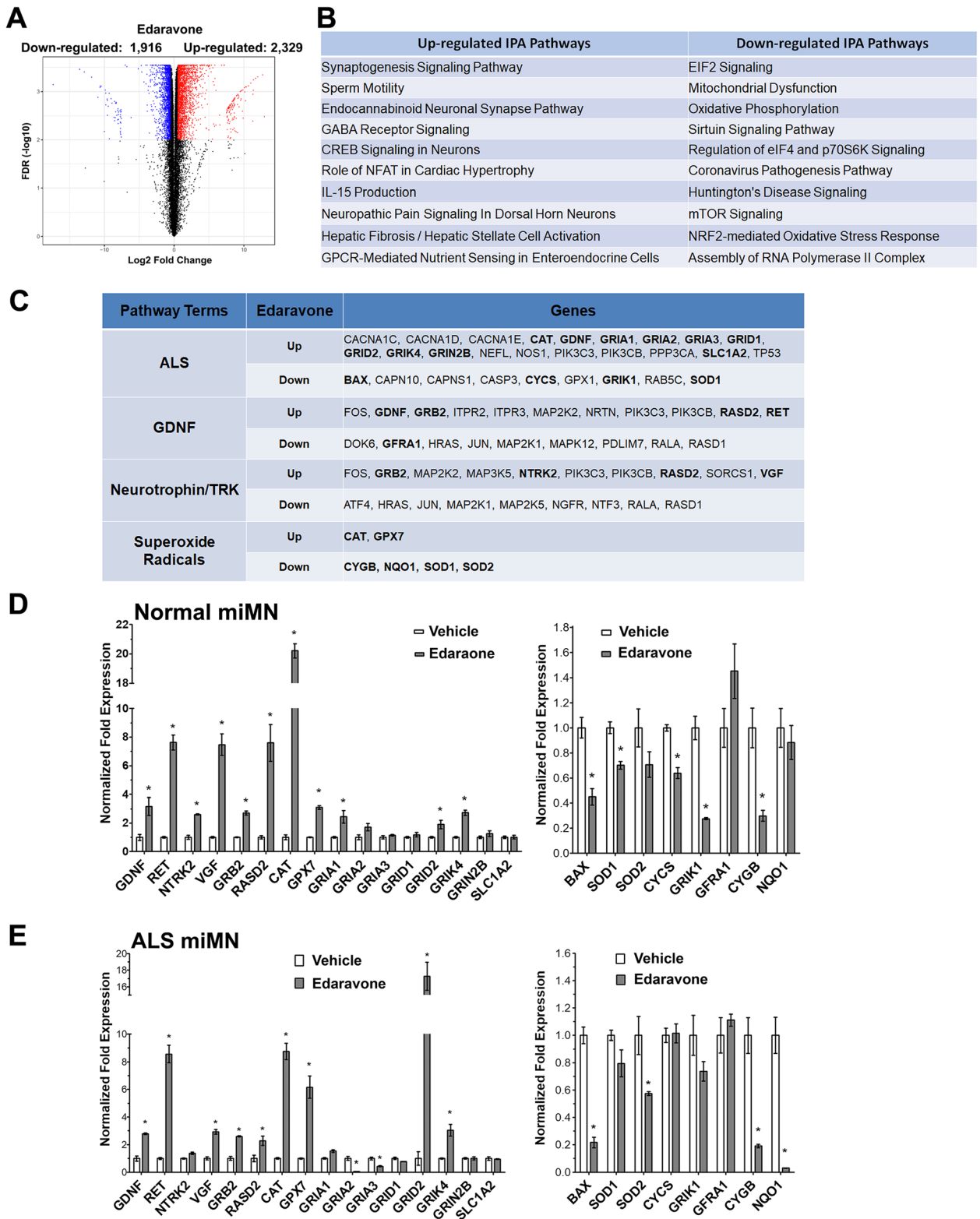


Fig. 6 (See legend on next page.)

(See figure on previous page.)

Fig. 6 Transcriptomic analysis of edaravone-induced responses in miMNs. **(A)** Volcano plot showing DE genes after edaravone treatment in control miMNs from iPSC1 (red and blue dots: genes with $FDR \leq 0.01$ and $\log_2(\text{fold change}) \geq 0.5$ or ≤ -0.5 , respectively). **(B)** Top ten IPA pathways enriched in up- and down-regulated genes. Pathways are ranked based on $-\log(p)$ value as calculated by the Fisher's exact test. **(C)** IPA pathways of interest and their associated genes. Genes selected for qPCR validation are highlighted. **(D and E)** Control and ALS miMNs were treated with edaravone (10 μM , 24 h) or DMSO as the vehicle control. Total cellular RNAs were subjected to qPCR analysis of up-regulated (left panel) and down-regulated (right panel) genes highlighted in C. Data represents Mean \pm SEM from 3 technical replicates (*: $p < 0.05$, t test with the Bonferroni correction for multiple comparisons)

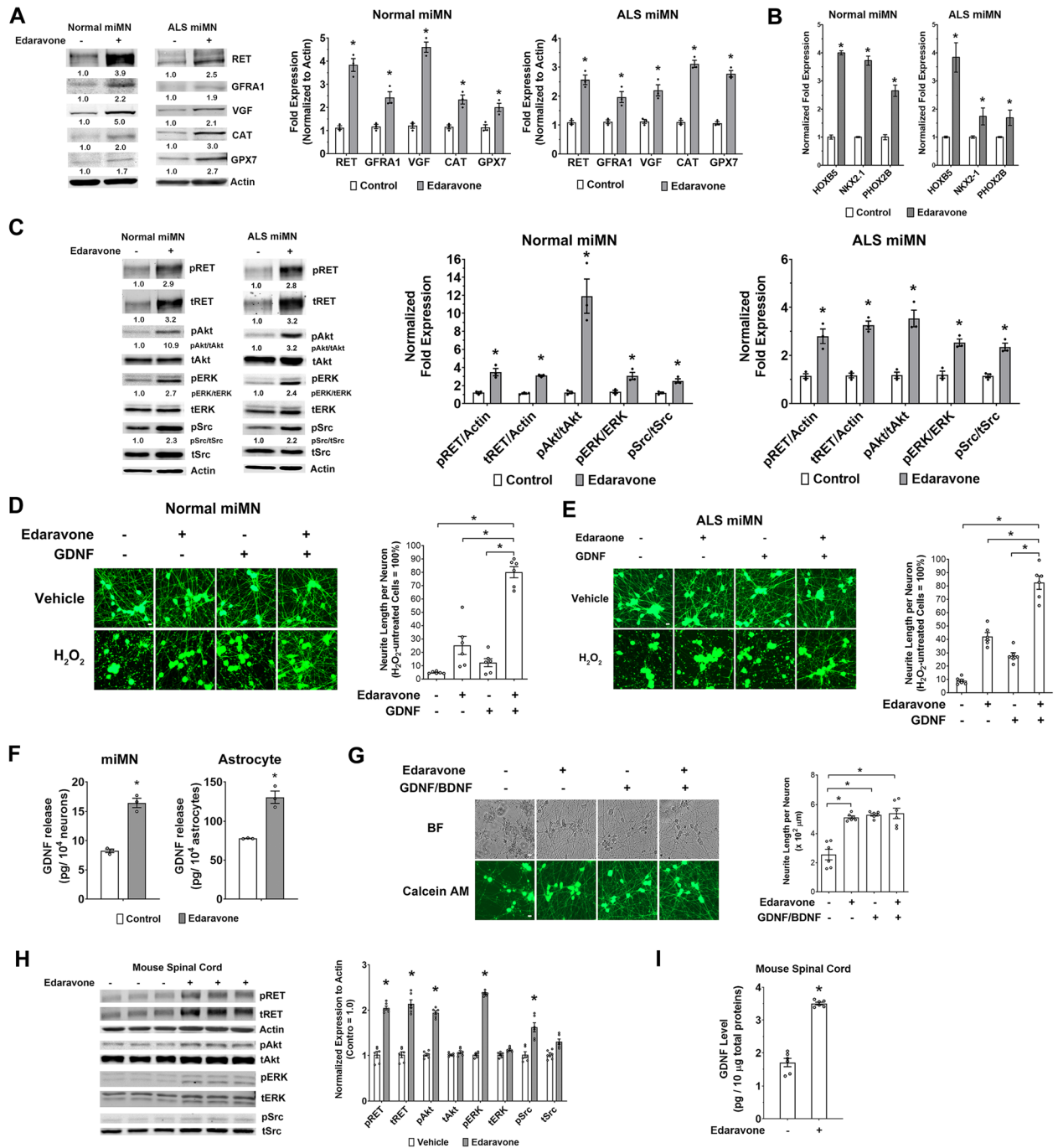


Fig. 7 (See legend on next page.)

(See figure on previous page.)

Fig. 7 Edaravone induces the GDNF/RET neurotrophic signaling pathway in miMNs and the mouse spinal cord. **(A and B)** Control and ALS miMNs were subjected to edaravone treatment (10 μ M, 24 h) or DMSO as the vehicle control. Total cellular proteins were subjected to western blotting of genes for validation (A, left) and quantification from three blotting results (B, right). Total cellular RNAs were subjected to qPCR analysis of three TF genes (B). (*: $p < 0.05$, t test with bonferroni correction for multiple comparisons). **(C)** Control and ALS miMNs with +/- edaravone treatment (10 μ M, 24 h) were subjected to western blotting (left) and quantification from three blotting results (right) to measure the levels of GDNF/RET signaling components (*: $p < 0.05$, t test with the Bonferroni correction for multiple comparisons). **(D and E)** Control and ALS miMNs (day 7 of differentiation) were pre-treated with +/- edaravone (10 μ M) and GDNF (1 ng/ml) for 16 h in neurotrophic factor-free medium and were treated with H₂O₂ (50 μ M) or PBS as the vehicle control for 24 h. Neurite length was quantified from 6 wells for each condition as technical replicates, after Calcium AM staining (Bar = 10 μ m). Edaravone+GDNF more effectively alleviated H₂O₂-induced neurite damage than single treatment. **(F)** Control miMNs and astrocytes differentiated from human NSCs were subjected to edaravone treatment (10 μ M) for 48 h. GDNF release in the culture medium was measured by ELISA with 3 technical replicates. **(G)** Control miMNs were cultured in neuron maturation medium with +/- edaravone and GDNF/BDNF for 15 days. Brightfield (BF) and Calcium AM staining images were shown (Bar = 10 μ m). Neurite length from 6 wells for each condition as technical replicates was compared among different conditions. **(H and I)** Mice ($n = 6$ for each group) received intraperitoneal injection of edaravone (15 mg/kg daily) or vehicle (saline) for 5 days. Total proteins from spinal cord tissues were analyzed by western blotting (G) Representative western blotting and quantification of protein fold expression normalized to Actin (*: $p < 0.05$, t test with the Bonferroni correction for multiple comparisons). Total proteins from spinal cord tissues from 6 mice were also analyzed by GDNF ELISA (H, *: $p < 0.01$, t test). Data represents Mean \pm SEM

miMNs showed higher levels of phosphorylated AKT, ERK and Src (Fig. 7C), three GDNF/RET-activated downstream kinases [43]. These results support that edaravone-induced RET expression leads to more active GDNF/RET signaling. We further asked if edaravone-induced RET expression leads to enhanced neuroprotection. We used a high dose of H₂O₂ treatment (50 μ M, 6 h) to induce severe neurite damage in control miMNs (Fig. 7D, about 95% neurite loss), which cannot be effectively protected by edaravone or GDNF alone (Fig. 7D, $p > 0.05$). In comparison, edaravone/GDNF combined treatment significantly inhibited H₂O₂-induced neurite damage (Fig. 7D, neurite loss: 37% compared to 95%, $p < 0.01$). Consistently, edaravone/GDNF combined treatment also protected ALS miMNs with the *SOD1*^{A4V} mutation (Fig. 7E, neurite loss: 18% compared to 91%, $p < 0.01$), more effectively than their single treatment (neurite loss: 58 and 72% for GDNF and edaravone single treatment, respectively).

RNA-seq and qPCR results both show that GDNF transcription is also induced by edaravone (Fig. 6). Consistently, we found that edaravone increased GDNF release from miMNs by about 100% (Fig. 7F, $p < 0.05$). In astrocytes differentiated from human fetal neural stem cells (NSCs), edaravone also increased GDNF release by 68% (Fig. 7F, $p < 0.05$). These results indicate that edaravone-induced GDNF release may lead to more active GDNF/RET neurotrophic signaling in autocrine and possibly also paracrine manners.

Since edaravone-induced RET expression and GDNF release may provide autocrine neurotrophic signaling to support neuron survival and maturation, we tested if edaravone can substitute the neurotrophic factors (GDNF and BDNF) used in miMN maturation medium. After 15 days of in-vitro maturation, we found that medium containing edaravone alone, GDNF/BDNF or both showed no difference in supporting normal

neuronal morphology and neurite outgrowth (Fig. 7G). miMNs from these three conditions also showed no difference in the expression of mature MN markers (e.g. ChAT and Synapsin 1) (data not shown). In contrast, miMNs survived poorly in medium without GDNF/BDNF or edaravone, and showed impaired neurite outgrowth (Fig. 7G). These results further support the neurotrophic function of edaravone-induced GDNF/RET signaling in miMNs.

Finally, we tested the in-vivo effect of edaravone on the GDNF/RET signaling in the spinal cord of mice with systemic edaravone treatment. Mice received intraperitoneal edaravone or vehicle treatment for 5 days. By analyzing total proteins from the entire spinal cord, we showed that edaravone treatment, compared to vehicle control, increased the levels of total and phosphorylated RET proteins by about 100%, and led to enhanced activation of RET downstream kinases (AKT, ERK and Src) (Fig. 7H). GDNF levels in mouse spinal cord tissues also showed more than 100% increase after edaravone treatment, as measured by ELISA (Fig. 7I). Overall, these results support edaravone's in-vivo efficacy to activate the GDNF/RET neurotrophic signaling.

Discussion

Synthetic mRNAs are highly efficient vehicles for delivering TF drivers to manipulate cell identity. Currently, mRNA-driven cell reprogramming kits are widely used for generating research and clinical grade human iPSCs [44]. Compared to cell reprogramming, mRNA-driven cell differentiation requires more optimization, including TF driver selection, TF sequence modification and mRNA delivery optimization, because a practical iPSC differentiation strategy requires at least 80% efficiency, compared to the 1–2% efficiency of cell reprogramming that is usually adequate [45, 46].

Traditional MN differentiation strategies for iPSCs use multiple compounds in a multi-step protocol and take 10–14 days to generate MNs with variable purity (commonly 50–70%) [15, 16]. Here, we established a 4-day one-step protocol to reproducibly generate miMNs with > 90% purity from iPSC lines from normal individuals and an ALS patient, as well as ESCs. We defined Ngn2 and Olig2 as a pair of TFs sufficient to drive highly efficient MN conversion. It is noteworthy to mention that Olig2 alone is not capable of inducing neuronal conversion. Compared to the Ngn2/Olig2 combination, a previous report used mRNAs coding five TFs including wild-type Ngn2 to generate human iPSC-derived MNs in 7 days [47]. Mazzoni et al. used three TFs (Ngn2, Isl1 and Lhx3) to differentiate mouse ESCs to MNs [18]. In the context of mRNA-based TF delivery, our results support that phosphosite modification in Ngn2 and Olig2 is a key element to increase the efficiency of mRNA-induced TF expression and MN conversion, as phosphosite modification leads to higher and long-lasting expression of Ngn2 and Olig2. SHH and DAPT were also found to promote MN induction, as MN lineage markers (HB9 and Islet1) were further induced by SHH/DAPT when N-SA/O-SA mRNAs were used. It is possible that mRNAs coding other TFs can be added to this protocol, which optimization may promote in-vitro survival and functional maturation of miMNs. Now, large-scale mRNA synthesis is easy and cost-effective. It is also noteworthy to mention that mRNAs used here were synthesized using nucleotides without pseudouridine and 5-methylcytidine modification, further reducing the cost of this method. In addition, miMNs after cryopreservation have a typical recovery rate of > 60%. Therefore, it is practical to establish a bank of iPSC-derived miMNs from familial and sporadic ALS patients, which platform will facilitate drug development for this devastating disease with complex biology and significant clinical heterogeneity [4]. We extensively characterized the identity of miMNs using well-defined MN markers as well as transcriptomic comparison with MNs differentiated by the traditional compound-based protocol [39]. Furthermore, we characterized their neuronal functions, using traditional patch-clamping recording as well as the MEA system suitable for high-throughput analysis. Moreover, high-content analysis of miMNs in the 1536-well format will allow ultra-high-throughput drug screening in an ALS MN bank, especially drug combination screening. It is noteworthy to mention that transcriptomic similarity between miMNs and compound-induced MNs at different maturation stages may be altered by batch effects in RNA-seq data. Thus, expression of mature neuron markers (e.g. Synapsin-1) and electrophysiological characterization may more reliably define the maturation status of miMNs.

Our following study of edaravone supports the applicability of miMNs in ALS drug development. Edaravone effectively protects miMNs from oxidative stress-associated neurotoxicity, consistent with its clinical efficacy shown in some ALS patients and proposed drug action as a free-radical scavenger to reduce oxidative stress. We also showed the neuroprotective function of edaravone in glutamate-induced neurotoxicity model using miMNs. Notably, the MEA analysis provides a more sensitive readout of neuronal damage as reflected by loss of neuronal activity. Since H₂O₂ is used at a much lower concentration without inducing rapid neurite damage, this MEA assay may more closely recapitulate a physiologically relevant condition of oxidative stress and neuronal dysfunction commonly seen during early ALS pathogenesis [4]. Compared to traditional screening platforms using neuron death or neurite damage as the readout, MEA analysis will help us identify potential compounds protecting neuronal functions of MNs and likely other neuron types involved in neurodegenerative disorders.

To the best of our knowledge, this study provides the first genome-wide map of edaravone-regulated signaling pathways in human MNs from iPSCs. This transcriptomic study will provide the foundation to explore mechanisms of action of edaravone and facilitate its clinical applications in various brain disorders. A significant discovery is edaravone-activated neurotrophic factor signaling pathways, especially GDNF/RET signaling. Edaravone induces the expression of GDNF receptors (e.g. RET) as well as GDNF release in miMNs, supporting by following results showing more active RET downstream kinases in edaravone-treated miMNs. Edaravone also induces GRB2 and RASD2, two key signaling molecules of the GDNF/RET signaling [43], which mechanism may further boost the activation of this neurotrophic signaling. GDNF is a potent neurotrophic factor promoting MN survival, and dysfunction of GDNF/RET signaling plays an essential role in ALS pathogenesis [48]. GDNF-based ALS therapies have been extensively studied in pre-clinical ALS models [49, 50], including recombinant GDNF proteins, GDNF gene therapy and GDNF-secreting cells. Some of these therapies have been advanced to human trials, such as GDNF-secreting neural progenitor cells (NCT02943850), but still need optimization. Our results demonstrate that the brain-penetrating ALS drug edaravone promotes RET expression and GDNF release in MNs, which mechanism likely mediates the neuroprotective effect of edaravone on miMNs. RET down-regulation has been found in MNs of the *SOD1*^{G93A} ALS mouse model. It is possible that ALS therapies simply supplying GDNF may be less effective than a therapeutic strategy that boosts both RET expression and GDNF supply. To support this

hypothesis, we showed that edaravone and GDNF combined treatment protected miMNs more effectively than their single treatment, which result justifies further pre-clinical and clinical studies of this combination strategy. Strikingly, edaravone can substitute all neurotrophic factors (GDNF/BDNF) in the miMN maturation medium to support in-vitro survival and maturation of miMNs, providing further evidence to support its neurotrophic function. To the best of our knowledge, this is the first report defining a small-molecule compound capable to replace neurotrophic factors essential for neuron culture. We also show that edaravone induces GDNF release in astrocytes, a major GDNF-secreting cell type in the brain, suggesting an edaravone-activated paracrine mechanism of neuroprotection. Moreover, edaravone shows consistent in-vivo efficacy to activate the GDNF/RET signaling in mouse spinal cord samples, justifying more studies in mouse models of ALS and other neurodegenerative diseases that may benefit from neuroprotection through GDNF/RET activation. Overall, our results provide a pharmacological strategy to activate the GDNF/RET neuroprotective signaling using a brain-penetrating drug with validated oral bioavailability.

The edaravone-induced neurotrophic protein VGF is a potential biomarker of ALS progression. VGF levels in the cerebral spinal fluid (CSF) and serum progressively decline in the *SOD1*^{G93A} ALS mouse model and ALS patients, and exogenous VGF is neuroprotective in MN cultures and ALS mouse models [51–53]. VGF is known to be induced by neurotrophic factors (e.g. BDNF and NGF) [54, 55] and also shows up-regulation after RET activation in neuroblastoma cells [56]. Thus, it is possible that edaravone induces VGF through the activation of GDNF/RET and/or BDNF/NTRK2 signaling. Elevated VGF levels in edaravone-treated miMNs may also function as a neuroprotective mechanism. Importantly, it is feasible to measure VGF-derived peptides in CSF and blood samples [51, 57]. Edaravone-induced VGF may be developed as a biomarker to predict drug response in clinical trials, such as the current ALS trial studying oral edaravone (NCT04165824).

Other edaravone-regulated genes with potential ALS association are also of interest, such as enzymes clearing superoxide radicals (e.g. CAT and GPX7), glutamate receptors (e.g. GRIA1, GRID2 and GRIK4) and pro-apoptotic proteins (e.g. BAX). More studies are required to understand how edaravone modulates various glutamate receptors that either positively or negatively regulate glutamate receptor-mediated excitotoxicity, a key pathogenetic mechanism in ALS [4, 58]. It is noteworthy to mention that our MEA analysis did not detect hyperexcitability in edaravone-treated miMNs.

Conclusion

We established a synthetic mRNA-driven strategy to efficiently generate iPSC-derived functional MNs applicable to high-throughput drug screening. We further define the neuroprotective effect of the ALS drug edaravone and reveal underlying molecular mechanisms, including the activation of GDNF/RET neurotrophic signaling. Methodology from this study will expand the applications of iPSC technology in MN research and therapeutic development for MN-associated neurodegenerative diseases. Novel molecular insights from this study will facilitate the development of edaravone-based therapies (e.g. edaravone/GDNF combination therapy) for ALS and likely other brain disorders associated with neurodegeneration (e.g. Parkinson's disease and stroke).

Abbreviations

iPSC: induced pluripotent stem cell; ESC: embryonic stem cell; PSC: pluripotent stem cell; MN: motor neuron; ALS: amyotrophic lateral sclerosis; miMN: mRNA-induced motor neuron; FDA: the Food and Drug Administration; TF: transcription factor

Supplementary Information

The online version contains supplementary material available at <https://doi.org/10.1186/s13024-021-00510-y>.

Additional file 1: Supplemental Fig. 1. (A) TUJ1+ miMNs from the iPSC1 line (day 7 of differentiation) show no detectable expression of the pluripotent stem cell marker (OCT4) and the oligodendrocyte lineage marker (O4). (B) TUJ1+ miMNs from the iPSC1 line were immunostained for the cholinergic neuron marker ChAT. Cytoplasmic signal of the ChAT protein was not detected in these miMNs at day 10 of differentiation, compared to cytoplasmic ChAT signal in more mature miMNs at day 20 of differentiation (Fig. 3A). (C) The iPSC line iPSC1 was differentiated by 3 daily transfections of NSA mRNA alone without using OSA mRNA and morphogenes. TUJ1+ NSA-induced neurons (day 30 of differentiation) express the glutamatergic neuron marker VGluT1, validating the VGluT1 antibody also used in Fig. 3E and supporting the requirement of Olig2 for MN induction. Cell nuclei were counterstained with DAPI (Bar = 20 μm). (D) Control miMNs from iPSC1 and iPSC3 iPSCs (referred to as control miMN 1 and 2, respectively) and ALS miMNs from iPSC2 and iPSC4 iPSCs (referred to as ALS miMN 1 and 2, respectively) were plated to the MEA plate at day 4 of differentiation. Their spontaneous spiking was recorded at indicated days in vitro (Fig. 4G). The number of active electrodes at indicated days in vitro did not show difference between control and ALS miMNs (3 technical replicates for each miMN line, linear regression with clustered data, ALS vs control miMNs). **Supplemental Fig. 2.** Reproducibility of two replicates from the transcriptomic analysis of miMNs with +/- edaravone treatment. (A) Heatmap clustering of RNA-Seq results from miMNs with +/- edaravone treatment (10 μM, 24 h, n = 2 for each group). Gene expression was calculated by reads per kilobase of transcript, per million mapped reads (RPKM). (B) A transcriptomic comparison between miMNs and differentiating cells at various days during compound-induced MN differentiation from human ESCs (Reference 39, GSE140747 from the GEO database). PCA plotting shows control (Con_1, Con_2) and edaravone-treated (Ed_1 and Ed_2) miMN samples (two technical replicates) more closely cluster with compound-induced MNs at day 8 and 9 of differentiation (three technical replicates labelled as D8_1/D8_2/D8_3 and D9_1/D9_2/D9_3, respectively). All samples are labelled as day at differentiation (D)_replicate number. All samples are grouped and colored based on the clustering result using model-based clustering of RNA-seq data.

Additional file 2: Table S1. Up-regulated genes in control miMNs after edaravone treatment.

Additional file 3: Table S2. Down-regulated genes in control miMNs after edaravone treatment.

Additional file 4: Table S3. Expression data of all genes detected by RNA-seq.

Additional file 5: Table S4. IPA signaling pathways enriched in edaravone up-regulated genes.

Additional file 6: Table S5. IPA signaling pathways enriched in edaravone down-regulated genes.

Additional file 7: Table S6. qPCR primers and antibodies.

Acknowledgements

We thank Dr. Ted Dawson and Dr. Valina Dawson for helping our studies of human pluripotent stem cells. We thank Dr. John Latterra for helpful discussion and editing the manuscript.

Authors' contributions

All authors read and approved the final manuscript. Q.L.: collection of data, data analysis and interpretation, manuscript writing, final approval of manuscript. Y.F.: collection of data, data analysis and interpretation, final approval of manuscript. Y.X.: collection of data, data analysis and interpretation, manuscript writing, final approval of manuscript. X.Z.: collection of data, data analysis and interpretation, manuscript writing, final approval of manuscript. Y.F.: collection of data, data analysis and interpretation, final approval of manuscript. G.G.: data analysis and interpretation, manuscript writing, final approval of manuscript. W.Z.: data analysis and interpretation, manuscript writing, final approval of manuscript. J.R.: provision of study material, final approval of manuscript. A.T.: provision of study material, final approval of manuscript. P.L.: provision of study material, final approval of manuscript. X.M.: provision of study material, final approval of manuscript. N.M.: conception and design, provision of study material, manuscript writing, final approval of manuscript. M.Y.: conception and design, financial support, collection of data, data analysis and interpretation, manuscript writing, final approval of manuscript. All authors read and approved the final manuscript.

Authors' information

None.

Funding

This work was supported by the Maryland Stem Cell Research Fund (M.Y. and N.M.).

Availability of data and materials

The RNA-Seq datasets reported here have been deposited to the Gene Expression Omnibus (GEO) Repository (GSE151997).

Declarations

Ethical approval and consent to participate

All experiments involving human stem cells have been approved by the Johns Hopkins Medicine Institutional Review Board. The animal protocol was approved by the Johns Hopkins School of Medicine Animal Care and Use Committee.

Consent for publication

Not applicable.

Competing interests

The authors declare that they have no competing interests.

Author details

¹Department of Endocrinology, Key Laboratory of Endocrinology, NHC, Peking Union Medical College Hospital, Chinese Academy of Medical Sciences & Peking Union Medical College, Beijing 100730, China. ²Hugo W. Moser Research Institute at Kennedy Krieger, 707 North Broadway, Baltimore, MD 21205, USA. ³Department of Physiology and Biophysics, Howard University, Washington, DC 20059, USA. ⁴Department of Biostatistics, Johns Hopkins Bloomberg School of Public Health, Baltimore, MD 21205, USA. ⁵Department of Psychiatry and Behavioral Sciences, Johns Hopkins University

School of Medicine, Baltimore, MD 21205, USA. ⁶Department of Neurology, Johns Hopkins University School of Medicine, Baltimore, MD 21205, USA.

Received: 11 June 2021 Accepted: 22 December 2021

Published online: 10 January 2022

References

- Qiang LFR, Abeliovich A. Remodeling neurodegeneration: somatic cell reprogramming-based models of adult neurological disorders. *Neuron*. 2013;78(6):957–69. <https://doi.org/10.1016/j.neuron.2013.06.002>.
- Ito DOH, Suzuki N. Accelerating progress in induced pluripotent stem cell research for neurological diseases. *Ann Neurol*. 2012;72(2):167–74. <https://doi.org/10.1002/ana.23596>.
- Jung YWHE, Kim KY, et al. Human induced pluripotent stem cells and neurodegenerative disease: prospects for novel therapies. *Curr Opin Neurol*. 2012;25(2):125–30. <https://doi.org/10.1097/WCO.0b013e3283518226>.
- Brown RH, Al-Chalabi A. Amyotrophic lateral sclerosis. *N Engl J Med*. 2017; 377(2):162–72. <https://doi.org/10.1056/NEJMra1603471>.
- Mejzini R, Flynn LL, Pitout IL, Fletcher S, Wilton SD, Akkari PA. ALS genetics, mechanisms, and therapeutics: where are we now? *Front Neurosci*. 2019;13: 1310. <https://doi.org/10.3389/fnins.2019.01310>.
- Hinchcliffe M, Smith A. Riluzole: real-world evidence supports significant extension of median survival times in patients with amyotrophic lateral sclerosis. *Degener Neurol Neuromuscul Dis*. 2017;7:61–70. <https://doi.org/10.2147/DNND.S135748>.
- Amyotrophic Lateral Sclerosis/Riluzole Study Group II ea. Dose-ranging study of riluzole in amyotrophic lateral sclerosis. *Lancet*. 1996;347(9013): 1425–31. [https://doi.org/10.1016/S0140-6736\(96\)91680-3](https://doi.org/10.1016/S0140-6736(96)91680-3).
- Miller RGMJ, Moore DH. Riluzole for amyotrophic lateral sclerosis (ALS) motor neuron disease (MND). *Cochrane Database Syst Rev*. 2012;3:1–34. <https://doi.org/10.1002/14651858.CD001447.pub3>.
- Abe K, Aoki M, et al. STE: safety and efficacy of edaravone in well defined patients with amyotrophic lateral sclerosis: a randomised, double-blind, placebo-controlled trial. *Lancet Neurol*. 2017;16(7):505–12. [https://doi.org/10.1016/S1474-4422\(17\)30115-1](https://doi.org/10.1016/S1474-4422(17)30115-1).
- Turnbull J. Reappraisal of an ALS trial: unaccounted procedural risk. *Lancet Neurol*. 2020;19(9):717–8. [https://doi.org/10.1016/S1474-4422\(20\)30265-9](https://doi.org/10.1016/S1474-4422(20)30265-9).
- Mora JS. Edaravone for treatment of early-stage ALS. *Lancet Neurol*. 2017; 16(10):772. [https://doi.org/10.1016/S1474-4422\(17\)30289-2](https://doi.org/10.1016/S1474-4422(17)30289-2).
- Akimoto M, Nakamura K. Writing group on behalf of the Edaravone ALSSG: Edaravone for treatment of early-stage ALS - Authors' reply. *Lancet Neurol*. 2017;16(10):772. [https://doi.org/10.1016/S1474-4422\(17\)30290-9](https://doi.org/10.1016/S1474-4422(17)30290-9).
- Takei K, Watanabe K, Yuki S, Akimoto M, Sakata T, Palumbo J. Edaravone and its clinical development for amyotrophic lateral sclerosis. *Amyotroph Lateral Scler Frontotemporal Degener*. 2017;18(sup1):5–10. <https://doi.org/10.1080/21678421.2017.1353101>.
- Rothstein JD. Edaravone: a new drug approved for ALS. *Cell*. 2017;171(4): 742. <https://doi.org/10.1016/j.cell.2017.10.011>.
- Maurly Y, Come J, Piskowski RA, Salah-Mohellibi N, Chevalyere V, Peschanski M, et al. Combinatorial analysis of developmental cues efficiently converts human pluripotent stem cells into multiple neuronal subtypes. *Nat Biotechnol*. 2015;33(1):89–96. <https://doi.org/10.1038/nbt.3049>.
- Du ZW, Chen H, Liu H, Lu J, Qian K, Huang CL, et al. Generation and expansion of highly pure motor neuron progenitors from human pluripotent stem cells. *Nat Commun*. 2015;6(1):6626. <https://doi.org/10.1038/ncomms7626>.
- Mizuguchi R, Sugimori M, Takebayashi H, Kosako H, Nagao M, Yoshida S, et al. Combinatorial roles of olig2 and neurogenin2 in the coordinated induction of pan-neuronal and subtype-specific properties of motoneurons. *Neuron*. 2001;31(5):757–71. [https://doi.org/10.1016/S0896-6273\(01\)00413-5](https://doi.org/10.1016/S0896-6273(01)00413-5).
- Mazzoni EO, Mahony S, Closser M, Morrison CA, Nedelec S, Williams DJ, et al. Synergistic binding of transcription factors to cell-specific enhancers programs motor neuron identity. *Nat Neurosci*. 2013;16(9):1219–27. <https://doi.org/10.1038/nn.3467>.
- Xue Y, Zhan X, Sun S, Karuppagounder SS, Xia S, Dawson VL, et al. Synthetic mRNAs drive highly efficient iPSC cell differentiation to dopaminergic neurons. *Stem Cells Transl Med*. 2019;8(2):112–23. <https://doi.org/10.1002/sctm.18-0036>.
- Sagal J, Zhan X, Xu J, Tilghman J, Karuppagounder SS, Chen L, et al. Proneural transcription factor Atoh1 drives highly efficient differentiation of

- human pluripotent stem cells into dopaminergic neurons. *Stem Cells Transl Med.* 2014;3(8):888–98. <https://doi.org/10.5966/sctm.2013-0213>.
21. De Filippis L, Lamorte G, Snyder EY, Malgaroli A, Vescovi AL. A novel, immortal, and multipotent human neural stem cell line generating functional neurons and oligodendrocytes. *Stem Cells.* 2007;25(9):2312–21. <https://doi.org/10.1634/stemcells.2007-0040>.
 22. Kim D, Paggi JM, Park C, Bennett C, Salzberg SL. Graph-based genome alignment and genotyping with HISAT2 and HISAT-genotype. *Nat Biotechnol.* 2019;37(8):907–15. <https://doi.org/10.1038/s41587-019-0201-4>.
 23. Pertea M, Pertea GM, Antonescu CM, Chang TC, Mendell JT, Salzberg SL. StringTie enables improved reconstruction of a transcriptome from RNA-seq reads. *Nat Biotechnol.* 2015;33(3):290–5. <https://doi.org/10.1038/nbt.3122>.
 24. Robinson MD, McCarthy DJ, Smyth GK. edgeR: a Bioconductor package for differential expression analysis of digital gene expression data. *Bioinformatics.* 2010;26(1):139–40. <https://doi.org/10.1093/bioinformatics/btp616>.
 25. Robinson MD, Oshlack A. A scaling normalization method for differential expression analysis of RNA-seq data. *Genome Biol.* 2010;11(3):R25. <https://doi.org/10.1186/gb-2010-11-3-r25>.
 26. Ohta Y, Nomura E, Shang J, Feng T, Huang Y, Liu X, et al. Enhanced oxidative stress and the treatment by edaravone in mice model of amyotrophic lateral sclerosis. *J Neurosci Res.* 2019;97(5):607–19. <https://doi.org/10.1002/jnr.24368>.
 27. Richner M, Jager SB, Siupka P, Vaegter CB. Hydraulic extrusion of the spinal cord and isolation of dorsal root ganglia in rodents. *J Vis Exp.* 2017. <https://doi.org/10.3791/55226>.
 28. Tanabe YRH, Jessell TM. Induction of motor neurons by sonic hedgehog is independent of floor plate differentiation. *Curr Biol.* 1995;5(6):651–8. [https://doi.org/10.1016/S0960-9822\(95\)00130-8](https://doi.org/10.1016/S0960-9822(95)00130-8).
 29. Roelink TQCH. The notch response inhibitor DAPT enhances neuronal differentiation in embryonic stem cell-derived embryoid bodies independently of sonic hedgehog signaling. *Dev Dyn.* 2007;236(3):886–92. <https://doi.org/10.1002/dvdy.21083>.
 30. Zhang Y, Pak C, Han Y, Ahlenius H, Zhang Z, Chanda S, et al. Rapid single-step induction of functional neurons from human pluripotent stem cells. *Neuron.* 2013;78(5):785–98. <https://doi.org/10.1016/j.neuron.2013.05.029>.
 31. Wang C, Ward ME, Chen R, Liu K, Tracy TE, Chen X, et al. Scalable production of iPSC-derived human neurons to identify tau-lowering compounds by high-content screening. *Stem Cell Reports.* 2017;9(4):1221–33. <https://doi.org/10.1016/j.stemcr.2017.08.019>.
 32. Akopian AN, Sivilotti L, Wood JN. A tetrodotoxin-resistant voltage-gated sodium channel expressed by sensory neurons. *Nature.* 1996;379(6562):257–62. <https://doi.org/10.1038/379257a0>.
 33. Khodakhah K, Melishchuk A, Armstrong CM. Killing K channels with TEA+. *Proc Natl Acad Sci U S A.* 1997;94(24):13335–8. <https://doi.org/10.1073/pnas.94.24.13335>.
 34. Sances S, Bruijn LI, Chandran S, Eggen K, Ho R, Klim JR, et al. Modeling ALS with motor neurons derived from human induced pluripotent stem cells. *Nat Neurosci.* 2016;19(4):542–53. <https://doi.org/10.1038/nn.4273>.
 35. Johnson MA, Weick JP, Pearce RA, Zhang SC. Functional neural development from human embryonic stem cells: accelerated synaptic activity via astrocyte coculture. *J Neurosci.* 2007;27(12):3069–77. <https://doi.org/10.1523/JNEUROSCI.4562-06.2007>.
 36. Takazawa T, Croft GF, Amoroso MW, Studer L, Wichterle H, Macdermott AB. Maturation of spinal motor neurons derived from human embryonic stem cells. *PLoS One.* 2012;7(7):e40154. <https://doi.org/10.1371/journal.pone.0040154>.
 37. Wainger BJ, Kiskinis E, Mellin C, Wiskow O, Han SS, Sandoe J, et al. Intrinsic membrane hyperexcitability of amyotrophic lateral sclerosis patient-derived motor neurons. *Cell Rep.* 2014;7(1):1–11. <https://doi.org/10.1016/j.celrep.2014.03.019>.
 38. Dell'Orco M, Milani P, Cereda C. Hydrogen peroxide-mediated induction of SOD1 gene transcription is independent from Nrf2 in a cellular model of neurodegeneration. *Biochim Biophys Acta.* 1859;2015(2):315–23. <https://doi.org/10.1016/j.bbaggm.2015.11.009>.
 39. Rayon T, Stamatakis D, Perez-Carrasco R, Garcia-Perez L, Barrington C, Melchionda M, et al. Species-specific pace of development is associated with differences in protein stability. *Science.* 2020;369(6510). <https://doi.org/10.1126/science.aba7667>.
 40. Mammen PP, Shelton JM, Ye Q, Kanatous SB, McGrath AJ, Richardson JA, et al. Cytochrome b5 is a stress-responsive hemoprotein expressed in the developing and adult brain. *J Histochem Cytochem.* 2006;54(12):1349–61. <https://doi.org/10.1369/jhc.6A7008.2006>.
 41. Zhu J, Garcia-Barcelo MM, Tam PK, Lui VC. HOXB5 cooperates with NKX2-1 in the transcription of human RET. *PLoS One.* 2011;6(6):e20815. <https://doi.org/10.1371/journal.pone.0020815>.
 42. Leon TY, Ngan ES, Poon HC, So MT, Lui VC, Tam PK, et al. Transcriptional regulation of RET by Nkx2-1, Phox2b, Sox10, and Pax3. *J Pediatr Surg.* 2009;44(10):1904–12. <https://doi.org/10.1016/j.jpedsurg.2008.11.055>.
 43. Airaksinen MS, Saarma M. The GDNF family: signalling, biological functions and therapeutic value. *Nat Rev Neurosci.* 2002;3(5):383–94. <https://doi.org/10.1038/nrn812>.
 44. Warren L, Manos PD, Ahfeldt T, Loh YH, Li H, Lau F, et al. Highly efficient reprogramming to pluripotency and directed differentiation of human cells with synthetic modified mRNA. *Cell Stem Cell.* 2010;7(5):618–30. <https://doi.org/10.1016/j.stem.2010.08.012>.
 45. Schlaeger TM, Dameron L, Brickler TR, Entwisle S, Chan K, Cianci A, et al. A comparison of non-integrating reprogramming methods. *Nat Biotechnol.* 2015;33(1):58–63. <https://doi.org/10.1038/nbt.3070>.
 46. Polegano MA, Eminli S, Beissert T, Herz S, Moon JJ, Goldmann J, et al. Efficient reprogramming of human fibroblasts and blood-derived endothelial progenitor cells using nonmodified RNA for reprogramming and immune evasion. *Hum Gene Ther.* 2015;26(11):751–66. <https://doi.org/10.1089/hum.2015.045>.
 47. Goparaju SK, Kohda K, Ibata K, Soma A, Nakatake Y, Akiyama T, et al. Rapid differentiation of human pluripotent stem cells into functional neurons by mRNAs encoding transcription factors. *Sci Rep.* 2017;7(1):42367. <https://doi.org/10.1038/srep42367>.
 48. Stanga S, Brambilla L, Kienlen-Campard P. A Role for GDNF and Soluble APP as Biomarkers of Amyotrophic Lateral Sclerosis Pathophysiology. *Front Neurol.* 2018;9:384.
 49. Wang LJ, Lu YY, Nakano I. Neuroprotective effects of glial cell line-derived neurotrophic factor mediated by an adeno-associated virus vector in a transgenic animal model of amyotrophic lateral sclerosis. *J Neurosci.* 2002;22(16):6920–8. <https://doi.org/10.1523/JNEUROSCI.22-16-06920.2002>.
 50. Thomsen GM, Avalos P, Ma AA, Alkaslasi M, Cho N, Wyss L, et al. Transplantation of neural progenitor cells expressing glial cell line-derived neurotrophic factor into the motor cortex as a strategy to treat amyotrophic lateral sclerosis. *Stem Cells.* 2018;36(7):1122–31. <https://doi.org/10.1002/stem.2825>.
 51. Zhao Z, Lange DJ, Ho L, Bonini S, Shao B, Salton SR, et al. Vgf is a novel biomarker associated with muscle weakness in amyotrophic lateral sclerosis (ALS), with a potential role in disease pathogenesis. *Int J Med Sci.* 2008;5:92–9. <https://doi.org/10.7150/ijms.5.92>.
 52. Noda Y, Motoyama S, Nakamura S, Shimazawa M, Hara H. Neuropeptide VGF-Derived Peptide LQEQ-19 has Neuroprotective Effects in an In Vitro Model of Amyotrophic Lateral Sclerosis. *Neurochem Res.* 2019;44(4):897–904. <https://doi.org/10.1007/s11064-019-02725-4>.
 53. Shimazawa M, Tanaka H, Hara H, et al. An inducer of VGF protects cells against ER stress-induced cell death and prolongs survival in the mutant SOD1 animal models of familial ALS. *PLoS One.* 2010;5(12):e15307. <https://doi.org/10.1371/journal.pone.0015307>.
 54. Mandolesi G, Gargano S, Pennuto M, Illi B, Molfetta R, Soucek L, et al. NGF-dependent and tissue-specific transcription of vgf is regulated by a CREB-p300 and bHLH factor interaction. *FEBS Lett.* 2002;510(1-2):50–6. [https://doi.org/10.1016/S0014-5793\(01\)03227-6](https://doi.org/10.1016/S0014-5793(01)03227-6).
 55. Alder J, Thakker-Varia S, Bangasser DA, Kuroiwa M, Plummer MR, Shors TJ, et al. Brain-derived neurotrophic factor-induced gene expression reveals novel actions of VGF in hippocampal synaptic plasticity. *J Neurosci.* 2003;23(34):10800–8. <https://doi.org/10.1523/JNEUROSCI.23-34-10800.2003>.
 56. Cerchia L, D'Alessio A, Amabile G, Duconge F, Pestourie C, Tavittian B, et al. An autocrine loop involving ret and glial cell-derived neurotrophic factor mediates retinoic acid-induced neuroblastoma cell differentiation. *Mol Cancer Res.* 2006;4(7):481–8. <https://doi.org/10.1158/1541-7786.MCR-06-0050>.
 57. Cocco C, Corda G, Lisci C, Noli B, Carta M, Brancia C, et al. VGF peptides as novel biomarkers in Parkinson's disease. *Cell Tissue Res.* 2020;379(1):93–107. <https://doi.org/10.1007/s00441-019-03128-1>.
 58. Leigh PN, Meldrum BS. Excitotoxicity in ALS. *Neurology.* 1996;47(Issue 6, Supplement 4):S221–7. https://doi.org/10.1212/WNL.47.6_Suppl_4.221S.

Publisher's Note

Springer Nature remains neutral with regard to jurisdictional claims in published maps and institutional affiliations.