SCIENTIFIC REPORTS

natureresearch

OPEN

Perfluorooctane sulfonate (PFOS) and perfluorooctanoate (PFOA) acutely affect human $\alpha_1\beta_2\gamma_{2L}$ GABA_A receptor and spontaneous neuronal network function *in vitro*

Anke M. Tukker, Lianne M. S. Bouwman, Regina G. D. M. van Kleef, Hester S. Hendriks, Juliette Legler & Remco H. S. Westerink^{*}

Concerns about the neurotoxic potential of polyfluoroalkyl substances (PFAS) such as perfluorooctane sulfonate (PFOS) and perfluorooctanoic acid (PFOA) increase, although their neurotoxic mechanisms of action remain debated. Considering the importance of the GABA_A receptor in neuronal function, we investigated acute effects of PFAS on this receptor and on spontaneous neuronal network activity. PFOS (Lowest Observed Effect Concentration (LOEC) 0.1 μ M) and PFOA (LOEC 1 μ M) inhibited the GABA-evoked current and acted as non-competitive human GABA_A receptor antagonists. Network activity of rat primary cortical cultures increased following exposure to PFOS (LOEC 100 μ M). However, exposure of networks of human induced pluripotent stem cell (hiPSC)-derived neurons decreased neuronal activity. The higher sensitivity of the $\alpha_1\beta_2\gamma_{2L}$ GABA_A receptor for PFAS as compared to neuronal networks suggests that PFAS have additional mechanisms of action, or that compensatory mechanisms are at play. Differences between rodent and hiPSC-derived neuronal networks highlight the importance of proper model composition. LOECs for PFAS on GABA_A receptor and neuronal activity reported here are within or below the range found in blood levels of occupationally exposed humans. For PFOS, LOECs are even within the range found in human serum and plasma of the general population, suggesting a clear neurotoxic risk.

Perfluorooctane sulfonate (PFOS) and perfluorooctanoic acid (PFOA) are well-known perfluoroalkyl substances (PFAS) consisting of an eight-carbon chain in which hydrogen atoms have been substituted with fluorine. Their combined hydrophobic, hydrophilic, oleophobic and lipophobic properties make them ideal industrial surfactants for the manufacturing of consumer products, including paint, stain repellents, fire-fighting foams, and non-stick cookware coatings. The strong carbon-fluorine bond renders PFOS and PFOA highly persistent and studies have shown their presence in the environment, wildlife and even human blood (for reviews see^{1,2}). By 2002, production of PFOS was phased out and production phase out of PFOA followed in 2006³. Recent studies indicate that these efforts may be responsible for a reduction in human blood levels in some areas, but the long half-lives of PFOS and PFOA result in slow elimination from environment and humans^{4,5}.

Research has demonstrated that the (developing) nervous system is one of the most sensitive targets for PFOS and PFOA. In mice and rats exposed pre- and/or neonatally to PFOS or PFOA, increased motor activity, decreased habituation and deficits in spatial learning and memory abilities have been observed⁶⁻¹¹. Developmental neuro-toxicity has also been observed in other species, including chicken¹² and zebrafish larvae^{13,14}. However, epide-miological studies have been inconclusive on the risks of PFAS exposure on neurodevelopment. Some studies indicate an association between prenatal PFOS and/or PFOA exposure and an increased risk on congenital cerebral palsy¹⁵, neuro-behavioural development¹⁶ and visual motor abilities¹⁷, whereas others reported no association between PFOS and PFOA exposure on psychomotor, cognitive and neurobehavioural development¹⁸, IQ levels¹⁹ or increased risk on ADHD and childhood autism²⁰. Notably, most early studies reported no or limited

Neurotoxicology Research Group, Toxicology Division, Institute for Risk Assessment Sciences (IRAS), Faculty of Veterinary Medicine, Utrecht University, P.O. Box 80.177, NL-3508, TD, Utrecht, The Netherlands. *email: R.Westerink@uu.nl

neurodevelopmental or neurological effects of PFAS exposure^{16,18,21-23}, although recent epidemiological studies indicate an association between PFAS exposure and child behaviour^{24,25}. It is important to keep in mind that exposure to PFAS in these studies is continuous and/or based on prolonged exposure paradigms.

The mechanism(s) underlying the observed neurotoxicity have been studied *in vitro* in neuronal cells. At (high) micromolar concentrations, PFOS and/or PFOA can affect a range of (pre)-synaptic processes, including reduction of cell viability^{26–29}, altered neuronal differentiation³⁰, increased formation of reactive oxygen species^{27,28,31–34}, and increased intracellular calcium (Ca²⁺) concentrations^{35,36}. In addition, PFOS and PFOA may also affect postsynaptic processes, including altered glutamate-activated currents and increase of potassium currents^{37,38}. By inducing influx of Ca²⁺ through voltage-dependent calcium channels, PFOS may exhibit acute excitotoxic effects on synaptic function and chronically inhibit synaptogenesis in rat brain hippocampal neurons³⁹. Although *in vivo* studies suggest the involvement of the cholinergic⁹ and glutamatergic system⁴⁰ in PFOS and PFOA mediated neurotoxicity, additional studies on postsynaptic receptors and channels by PFOS or PFOA are lacking.

Previous studies have shown agonistic effects on the postsynaptic human GABA_A receptor by several persistent organic pollutants, including polychlorinated biphenyls and brominated flame retardants^{41–44}. The GABA_A receptor is the main inhibitory neurotransmitter receptor in the central nervous system and is critical for brain development⁴⁵, long-term potentiation and synaptic plasticity⁴⁶, but also for functional neuronal signal transduction. Because the GABA_A receptor is the most important inhibitory receptor in the central nervous system, any effect of PFAS on this receptor may in turn affect neuronal signal transduction and spontaneous neuronal network activity.

In the present study we therefore used *Xenopus* oocytes that express the human $\alpha_1\beta_2\gamma_{2L}$ GABA_A receptor, which is the most abundant subunit combination⁴⁷, to investigate the acute effects of PFOS and PFOA using the two-electrode voltage-clamp technique. Additionally, we investigated whether the effects of PFOS and PFOA on the $\alpha_1\beta_2\gamma_{2L}$ GABA_A receptor are reflected in the level of spontaneous neuronal network activity using micro-electrode array (MEA) recordings in rat primary cortical neuronal networks, the current gold standard for MEA assays. These experiments were followed by measurements in human iPSC-derived neuronal models to investigate whether they were similarly affected by PFAS exposure as the rodent model.

Results

Antagonistic effects of PFAS on human GABA_A receptor. Oocytes expressing the human α_1 , β_2 , γ_{2L} subunits were superfused with various concentrations of GABA to derive a concentration-response curve with an EC₂₀, EC₅₀ and Hill slope of 36 μ M, 103 μ M and 1.3, respectively (see Fig. S1 and in line with previous publications⁴⁸). The expression and functionality of the γ_{2L} subunit was demonstrated by co-exposing GABA-responsive oocytes to a GABA-diazepam co-exposure (see Fig. S1).

To investigate whether PFOS and PFOA are able to activate the human GABA_A receptor, GABA-responsive oocytes were superfused with saline containing 0.1–100 μ M PFOS or PFOA. Neither PFOS nor PFOA induced any ion current, demonstrating that PFOS and PFOA by themselves are not agonists of the GABA_A receptor (data not shown). To determine possible partial agonistic or antagonistic effects, PFOS and PFOA were co-applied with a low concentration of GABA (~EC₂₀) and the lowest observed effect concentrations (LOEC; defined as the lowest test concentration that is significantly different from the control) of PFOS and PFOA were calculated. At this low effective GABA concentration, both PFOS (LOEC 0.1 μ M, IC₅₀ 0.28 μ M, CI 0.22–0.34 μ M, see Fig. 1a,e and Table 1; F(1,21) = 10.2, *p* = 0.004) and PFOA (LOEC 10 μ M, IC₅₀ 0.28 μ M, CI 19.28–24.21 μ M, see Fig. 1b,f and Table 1; F(1,22) = 256, *p* = <0.0001) concentration-dependently inhibited the GABA-evoked ion current. These data demonstrate that both PFOS and PFOA exert antagonistic effects on the GABA_A receptor. Notably, as shown in Fig. 1a, the PFOS-induced inhibition of the GABA-evoked ion current is not readily reversible. In contrast, the PFOA-induced inhibition is reversed within seconds (Fig. 1b).

Next, PFOS and PFOA were co-applied with a high concentration GABA (1 mM, EC₁₀₀). PFOS induced a poorly reversible, concentration-dependent inhibition of the GABA-evoked ion current (LOEC 0.1 μ M, IC₅₀ 0.18 μ M, CI 0.15–0.22 μ M; F(1,20) = 10.4, *p* = 0.004; Fig. 1c,e and Table 1), whereas the concentration-dependent PFOA-induced inhibition is rapidly reversed (LOEC 1 μ M, IC₅₀ 6.7 μ M, CI 5.96–7.59 μ M; F(1,22) = 66.2, *p* = <0.0001; Fig. 1d,f and Table 1).

The combined data indicate that PFOS and PFOA are potent, non-competitive GABA_A receptor antagonists. As such, PFOS and PFOA may antagonize the inhibitory function of the CNS. To determine whether the effects of these PFAS also modulate neuronal network function, effects of PFOS and PFOA on spontaneous neuronal network activity were assessed.

Effects of PFAS on rat cortical network activity. Recently, non-invasive measurements of neuronal activity using MEA recordings have shown that rat primary cortical cultures develop spontaneous network activity⁴⁹, which can be modulated by a diverse range of chemicals^{50–53}. Currently, rat primary cortical neurons are the gold standard for MEA recordings. At DIV9, network bursts throughout the well can be observed. Therefore, we exposed rat primary cortical cultures at DIV9–11 to PFOS or PFOA to assess effects on spike, burst and network burst related parameters.

Acute exposure to PFOS or PFOA affected mean spike rate (MSR; Fig. 2A), with statistically significant differences between the concentrations (Welch's F(4, 44.7) = 9.76, p = <0.0001 for PFOS and Welch's F(4, 46.5) = 3.5, p = 0.01 for PFOA). Following exposure to 100 µM PFOS, a marked increase in MSR could be observed (175 ± 13.1%; p = <0.0001). PFOS also increased the mean burst rate (MBR; Fig. 2B), with a statistically significant difference between concentrations (Welch's F(4, 46.0) = 7.2, p = 0.0001). However, this increase was only significant compared to control following exposure to 100 µM PFOS (175 ± 16.7%; p = <0.0001). PFOS did not affect mean network burst rate (MNBR; Fig. 2C). PFOA only induced minor effects on spike and (network) burst



Figure 1. Antagonistic effects of PFOS and PFOA on the human GABA_A receptor. Example recordings of inhibition of GABA-evoked ion currents by co-application of PFOS (1 μ M, **a**, **c**) or PFOA (10 μ M, **b**, **d**) with GABA (at low and high effective GABA concentration, respectively). PFOS-induced inhibition is poorly reversible (**a**, **c**), whereas PFOA-induced inhibition is reversed within seconds (**b**, **d**). Scale bar applies to all traces. Concentration-response curves show the concentration-dependent inhibition of GABA-evoked responses by PFOS (**e**) and PFOA (**f**) on the human GAB_A receptor with at low (**e**,**f**; solid line) and high (**e**,**f**; dashed line) effective GABA concentrations. Inhibition is presented as percentage of the GABA-evoked response (mean ± SEM, n = 3-4 oocytes per concentration from N = 1-2 batches).

rate. PFOA increased MNBR at 10μ M, but this increase was not significant. Neither PFOS nor PFOA affected burst duration (Fig. 2D), although there is a significant difference between groups for PFOA (Welch's F(4, 45.8) = 3.81, p = 0.009). Both PFAS did not affect network burst duration (Fig. 2E), but for PFOA the one-way Welch ANOVA was significant (Welch's F(4, 45.9.) = 3.67, p = 0.011). Effects observed at the highest test concentration are not a result of cytotoxicity (see Supplemental Data Fig. S2).

Overall, PFOS has a more profound effect on spontaneous neuronal network activity than PFOA. However, these data indicate that PFOS and PFOA induce only mild hyperexcitation in rat primary cortical cultures, which was unexpected considering the strong inhibition of human GABA_A receptors. This apparent discrepancy may be the result of inter-species differences. To investigate whether this is the case, human induced pluripotent stem cell (hiPSC)-derived neurons can be used. It is known that there is variability between different hiPSC-derived neuronal cell lines⁵⁴. This could be the result of the generation protocol or the maturation stage in which the neurons were frozen. To take these potential differences into account, two commercially available hiPSC-derived neuronal cell models were exposed to PFOS or PFOA: the iCell Glutaneuron – iCell Astrocytes co-culture and the SynFire iNS co-culture.

		Low effective GAB	A concentration	High effective GABA concentration		
	Concentration (µM)	% Inhibition	p-value	% Inhibition	p-value	
PFOS	0.01	4.8 ± 4.8 (4)	n.s	7.6±3.3 (4)	n.s	
	0.1	15 ± 1.8 (4)	0.002	33±4.6 (4)	< 0.0001	
	1	90±1.5(3)	< 0.0001	89±0.9(3)	< 0.0001	
	10	98±1.2(4)	< 0.0001	100±0(3)	< 0.0001	
	100	100 ± 0 (4)	< 0.0001	100±0(4)	< 0.0001	
	0.01	0±0(4)	n.s.	0.4±0.5 (4)	n.s	
PFOA	0.1	0±0(4)	n.s.	0±0(3)	n.s	
	1	6 ± 0.4 (4)	n.s	16±0.4 (4)	< 0.0001	
	10	31±1.9 (4)	< 0.0001	58±1.5(4)	< 0.0001	
	100	83±2.7 (4)	< 0.0001	93±2.7 (4)	< 0.0001	

Table 1. Inhibition of GABA-evoked ion current by PFOS or PFOA. Mean percentage of inhibition of the current (\pm SEM (*n*)) induced during exposure to PFOS or PFOA. Inhibitions were compared to control and *p*-values are given; n.s.: non-significant.

Effects of PFAS on spontaneous hiPSC network activity. Recent work in our laboratory has shown that hiPSC-derived neuronal co-cultures develop spontaneous activity and (network) bursting behaviour⁵⁵. Acute exposure to PFOS (Welch's F(4, 34.9) = 278, $p = \langle 0.0001 \rangle$ and PFOA (Welch's F(4, 32.9) = 13, $p = \langle 0.0001 \rangle$ affected MSR in the iCell Glutaneuron - iCell Astrocytes co-culture (Fig. 3A)). PFOS significantly decreased the MSR at $0.1 \,\mu$ M (91.8 ± 3.18%; *p* = 0.04) and 100 μ M (8.2 ± 3.71%; *p* = <0.0001). For PFOA, MSR was significantly decreased at 1 μ M (88.9 \pm 2.64; p = 0.02), 10 μ M (87.4 \pm 2.5%; p = 0.02) and 100 μ M (59.7 \pm 6.49; p < 0.0001) as compared to the control. Bursting was significantly affected following exposure to either PFAS (Welch's F(4, 34.9) = 36.0, p < 0.0001 for PFOS and Welch's F(4, 36.7) = 6.67, p = 0.0004; Fig. 3B). PFOS exposure significantly decreased MBR at $100 \,\mu\text{M}$ (9.43 \pm 4.2%; p = <0.0001). For PFOA, MBR significantly decreased at $10 \,\mu\text{M}$ (59.8 $\pm 10.7\%$; p = 0.023) and $100 \,\mu\text{M}$ (49.5 $\pm 8.73\%$; p = 0.0035). MNBR was decreased following exposure to PFOS (Welch's F(4, 30.1) = 51.0, $p = \langle 0.0001 \rangle$ and PFOA (Welch's F(4, 35.4) = 15.6, $p = \langle 0.0001 \rangle$; Fig. 3C). The decrease was significant following exposure to PFOS at $0.1 \,\mu\text{M}$ ($54.8 \pm 9.1\%$; p = 0.002) and $100 \,\mu\text{M}$ $(2.58 \pm 1.9\%; p < 0.0001)$ and for PFOA at 1 μ M ($65.4 \pm 10.1\%; p = 0.01$), 10 μ M ($44.7 \pm 9.03; p = 0.0005$) and $100 \,\mu\text{M}$ (22.2 ± 6.11; p < 0.0001). Exposure of the iCell Glutaneuron – iCell Astrocytes co-culture to either of the PFAS increased burst duration (Welch's F(4, 24.0) = 6.74, p = 0.0009 for PFOS and Welch's F(4, 31.6) = 6.71, p = 0.0005 for PFOA; Fig. 3D). The increase in burst duration following exposure to PFOS was significant at $10 \,\mu\text{M}$ (206 ± 39.3; p = 0.002) and $100 \,\mu\text{M}$ (245 ± 41.3%; p = 0.0002). Following exposure to PFOA, the increase was significant at 1 μ M (194 \pm 18.9%; p = <0.0001) and 100 μ M (174 \pm 30.6; p = 0.01). There was no significant effect on network burst duration (Fig. 3E), despite an apparent increase following exposure to 10 µM PFOS. This change in activity pattern is further illustrated by spike raster plots that show the decrease in spiking and (network) bursting activity combined with prolonged (network) burst duration following exposure to 10 µM PFOS (Fig. 4A) or 100 µM PFOA (Fig. 4B).

Overall, effects are more pronounced following exposure to PFOS. However, PFOS and PFOA do not induce hyperexcitation in the iCell Glutaneuron – iCell Astrocytes co-culture model. Since the chemical sensitivity can differ between different hiPSC-derived neuronal cultures⁵⁶, an additional hiPSC-derived cell model was tested, the SynFire iNS co-culture.

The SynFire iNS co-culture contains a higher ratio of astrocytes compared to the iCell Glutaneuron – iCell Astrocytes co-culture and although the ratio glutamatergic to GABAergic neurons is comparable, the total cell number is much higher thus the number of glutamatergic neurons is also higher in this culture⁵⁶. Therefore, the SynFire iNS co-culture is more active and excitable. Exposure of SynFire iNS co-culture to PFOS affected the MSR (Welch's F(4, 16.9) = 3.18, p = 0.04 for PFOS) with a significant decrease at 100 µM (69.1 ± 12; p = 0.04; Fig. 5A) as compared to control. Exposure to PFOA decreased the MSR at 100 µM, but this decrease was not significant. MBR was only significantly decreased by PFOS (Welch's F(4, 18.2) = 5.24, p = 0.006) at 100 µM (58.3 ± 8.13; p = 0.04; Fig. 5B). Neither PFOS nor PFOA significantly disturbed MNBR (Fig. 5C). Burst duration and network burst duration were not significantly affected, although an increase could be observed following exposure to PFOS 100 µM (Fig. 5D,E).

Comparison of PFAS effects in different cell models. LOECs for the activity metrics MSR, MBR and MNBR differ between the three different cell models (Table 2). The iCell Glutaneuron – iCell Astrocytes co-culture is the only model for which LOECs could be defined for PFOS and PFOA on all three metrics. This hiPSC-derived neuronal model is also the most sensitive of the three. The rat primary cortical neurons and SynFire iNS co-culture model show comparable LOECs, but respond differently to PFOS and PFOA exposure.

To further illustrate the effect of PFOS and PFOA on the rat primary cortical neurons and the two hiPSC-derived neuronal models, we created a heatmap including 19 parameters that are descriptive for neuronal (network) activity (Fig. 6; Table 3). Although many changes (and some concentration-dependent changes) are visible in the heatmap and in Fig. 3, it should be noted that many of these changes do not reach statistical significance. Nevertheless, it is clear that the rat cortex is rather insensitive, though an increased activity can be observed following exposure to 100 µM. The iCell Glutaneuron – iCell Astrocytes co-culture model is the most sensitive



Figure 2. Modulation of spontaneous neuronal activity of rat primary cortical neurons exposed to PFOS (black) or PFOA (grey) following acute exposure. Effects on mean spike rate (MSR; **A**), mean burst rate (MBR; **B**), mean network burst rate (MNBR; **C**), burst duration (**D**) and network burst duration (**E**) are depicted as average in % change of control (solvent control set to 100%; dashed line) \pm SEM from n = 18–20 wells and N = 4 plates. *Indicates p < 0.05.



Figure 3. Modulation of spontaneous neuronal activity of iCell Glutaneuron – iCell Astrocytes co-culture exposed to PFOS (black) or PFOA (grey) following acute exposure. Effects on mean spike rate (MSR; **A**), mean burst rate (MBR; **B**), mean network burst rate (MNBR; **C**), burst duration (**D**) and network burst duration (**E**) are depicted as average in % change of control (solvent control set to 100%; dashed line) \pm SEM from n = 3-17 wells and N = 3-5 plates. *Indicates p < 0.05.

model. This can be seen from the profound decrease in MSR, MBR and MNBR and the increase in burst duration. The SynFire iNS co-culture is just as (in)sensitive as the rat primary cortical culture (Figs. 2, 5 and 6), though with an increase, in contrast to a decrease, in activity following exposure to $100\,\mu$ M. It is also apparent that PFOS has a stronger and more excitatory effect than PFOA. In all three models, synchronicity parameters are least affected. Overall, this highlights that there are differences between human and rodent models. It also indicates that there



Figure 4. Spike raster plots illustrating the pattern of activity of a representative well of the iCell Glutaneuron – iCell Astrocytes co-culture before exposure (left) and the same well following exposure (right) to PFOS $10 \,\mu$ M (**A**) or PFOA $100 \,\mu$ M (**B**). Each row depicts one electrode in a well and each tick mark represents one spike in a 50 s interval. Spikes are depicted in black and bursts are depicted in blue.

are differences between the human model systems. Overall the model system used can to a large degree determine the type and degree of the effect.

Discussion

PFOS and PFOA are omnipresent in the environment. In vivo and in vitro research have indicated the neurotoxic potential of both compounds⁵⁷. However, the underlying mechanisms of action are still debated. In the present study we show for the first time that PFOS and PFOA exhibit concentration-dependent antagonistic effects on the human $\alpha_1\beta_2\gamma_{2L}$ GABA_A receptor at sub-micromolar concentrations (Fig. 1, Table 1). It is important to note that we tested only the most abundant $\alpha_1\beta_2\gamma_{2L}$ receptor subunit combination. As many other GABA_A receptor subtypes are present in the nervous system, PFAS may have different effects on these different receptor subtypes. The GABA-evoked current is rapidly restored once exposure to PFOA stops, but reversibility of the effect caused by PFOS is slow. This indicates that PFOA most likely has a lower receptor affinity than PFOS, which could be caused by the differences in head group. The reversibility could explain why at a high effective GABA concentration the LOEC of PFOA (1 µM) is tenfold higher than the LOEC of PFOS (0.1 µM). The observed inhibition of the GABA-evoked current occurs independently of the effective GABA concentration, implying that both PFAS are non-competitive GABA antagonists. This also indicates that neither PFOS nor PFOA binds to the GABA binding site located at the interface between the α and β subunit, but are more likely to bind to modulatory binding sites. A variety of compounds, including convulsants and barbiturates, target these modulatory binding sites⁵⁸. However, when positive or negative allosteric modulators bind to modulatory binding sites, the effect of the compound usually decreases with an increasing concentration of GABA. This is in contrast with our results. The agonistic effects of PFOS and PFOA could therefore be the result of a pore block through insertion of (a part of) the molecule into the receptor pore. This pore block would then prevent influx of chloride ions following binding of GABA to the GABA binding site. Full elucidation of the precise mode of action of PFAS on $\alpha_1\beta_2\gamma_{2L}$ GABA_A receptor function would require extensive mechanistic research, including measuring voltage-current curves in presence and absence of PFAS. Such experiments may elucidate whether PFOS and/ or PFOA cause a voltage-dependent pore block, which is a likely mechanism of action given our results. However, regardless the exact molecular mechanism, the antagonistic effects of PFOS and PFOA on the human GABA_A receptor can be hypothesized to cause hyperexcitation in neuronal networks.

This hypothesis is further supported by evidence that PFOS induces the influx of extracellular calcium³⁹. However, when primary rat cortical neurons are exposed to PFAS, excitation is observed only in spike and burst activity at 100 μ M PFOS, whereas no significant excitatory effects of PFOA were found (Figs. 2 and 6). As there were no effects on primary rat cortical neurons at low or submicromolar levels of PFAS, as observed for human



Figure 5. Modulation of spontaneous neuronal activity of SynFire iNS co-culture with PFOS (black) or PFOA (grey). Effects on mean spike rate (MSR; **A**), mean burst rate (MBR; **B**), mean network burst rate (MNBR; **C**), burst duration (**D**) and network burst duration (**E**) are depicted as average in % change of control (solvent control set to 100%; dashed line) \pm SEM from n = 7-9 wells and N = 2 plates. *Indicates p < 0.05.

	MSR		MBR		MNBR	
	PFOS	PFOA	PFOS	PFOA	PFOS	PFOA
Rat primary cortical neurons	100 µM	_	100 µM	—	_	—
iCell Glutaneuron – iCell Astrocytes co-culture	0.1 µM	1 µM	100 µM	10 µM	0.1 µM	1 µM
SynFire iNS co-culture	100 µM	—	100 µM	—	—	—

Table 2. LOECs of PFOS and PFOA for the different activity parameters (mean spike rate (MSR), mean burst rate (MBR) and mean network burst rate (MNBR)) on the different cell models. – Indicates no LOEC could be defined.

GABA_A receptor, two different hiPSC-derived neuronal co-cultures were exposed to PFOS and PFOA to exclude inter-species differences. In contrast to the expected hyperexcitation, neither of the hiPSC-derived neuronal models showed hyperexcitation following acute exposure to PFOS or PFOA. In the iCell Glutaneurons - iCell Astrocytes co-culture, a strong inhibition of activity in combination with a prolonged burst duration was found following exposure to PFOS or PFOA (Figs. 3, 4 and 6). This model showed spike and network burst LOECs for PFOS (0.1 μ M) and PFOA (1 μ M) that were comparable to the human GABA_A receptor studies. In contrast, in the SynFire iNS co-culture showed an apparent lack of sensitivity comparable to the rat primary cortical culture (Figs. 5, 6). While the human iPSC models thus differ in sensitivity, they both show inhibition of neuronal activity in contrast to the rat primary culture. This hints towards a specific species-difference. However, giving the inhibitory effect on the human $\alpha_1\beta_2\gamma_{2L}$ GABA_A receptor, these findings also imply that PFAS may have additional targets that mask the inhibition of the GABA_A receptor and/or that compensatory mechanisms are at play in neuronal networks to dampen the loss of GABA_A receptor function. For example, PFAS could act on different neurotransmitter receptors, such as glutamate receptors, thereby masking the inhibition of the GABA_A receptor. This idea is supported by research that showed that PFOS induces NMDA receptor mediated excitotoxicity⁵⁹. Toxicity results of a high influx of extracellular Ca^{2+} , which in its turn over-activates the NMDA receptor. Although this study supports the argument that PFAS hit multiple targets, it should be noted that these results were obtained following 24 h exposure, as compared to our acute exposure study. Alternatively, the hyperexcitability effects of PFOS and PFOA may become apparent only after chronic exposure. These effects may then be the result of changes in for example the AMPA receptor composition or expression, as it has been previously shown that prolonged exposure of rat primary cortical neurons to 1 µM PFOS lowers the expression of the AMPA receptor subunit GluR2, resulting in increased intracellular calcium concentrations and enhanced glutamate sensitivity⁶⁰. It is also possible that the differences between the rodent and human neuronal model are due to an immature phenotype of the hiPSC-models. However, the used models have been shown to have a mature phenotype with hyperpolarizing GABA-responses and expression of mature transporters like KCC2 and NKCC^{61,62}.



Figure 6. Heatmap of the effects of PFOS (light green) and PFOA (darker green) on selected metric parameters on rat primary cortical neurons (top; grey), iCell Glutaneuron – iCell Astrocytes (middle; black) and SynFire iNS co-culture (bottom; green). Colour scaling is based on the magnitude change in percentages relative to vehicle control based on n = 18-20 wells and N = 4 plates for rat primary cortical neurons; n = 3-17 wells and N = 3-5 plates for iCell Glutaneuron – iCell Astrocytes; n = 6-9 wells N = 2 plates for SynFire iNS co-culture.

Compared to the two hiPCs-derived neuronal models, the rat primary cortical culture contains the highest ratio of astrocytes, about 45%^{63,64}. Astrocytes increase the level of network activity^{55,65,66} and modulate responses to toxicological insults⁶⁷. It has been shown that astrocytes can also provide protection against adverse effects of PFOS on rat hippocampal neurons⁶⁸. Therefore, the lack of an excitatory effect at lower concentrations could be due to the protective effect of the astrocytes present. When astrocytes are exposed for 24 h to concentrations higher than 50 µM PFOS, astrotoxicity starts to occur, thereby increasing the risk on excitotoxicity⁶⁸. Although exposure duration in our study was shorter and no cytotoxicity was observed, the increase in spike and burst activity following acute exposure of rat primary cortical neurons to a high PFOS concentration could be indicative of the onset of excitotoxicity. Another potential explanation for the inhibition of activity in the hiPSC-derived neuronal cultures could reside in the PFOS-induced shift in the activation curves of voltage-gated sodium channels towards hyperpolarization^{37,69}, making it harder for the cell to reach an action potential. Future research to elucidate the possible counteracting mechanisms of action observed in our network study of PFOS and PFOA could focus on patch-clamp studies in neuronal networks following both acute and chronic exposure.

For PFOS and PFOA human serum to plasma ratios are $1:1^{70}$. Reported plasma or serum concentrations for PFOA fall for the general population in the range between $0.001-0.2 \ \mu M^{71-73}$. In the occupationally exposed population, levels as high as $2.5 \ \mu M$ are found⁷⁰. Thus, LOECs reported on GABA_A receptor function, spike rate and network burst activity ($1 \ \mu M$) are below this level. Importantly, PFOA can cross the blood-brain barrier⁷². Therefore, PFOA could be a risk for the occupationally exposed population. PFOS can cross the blood-brain barrier⁷², and or serum concentrations for PFOS fall in the range between $0.002-1.3 \ \mu M$ in the general population⁷¹⁻⁷³ and in occupationally exposed workers the levels are even higher, ranging from $0.27 \ to 10 \ \mu M^{70,71}$. LOECs on GABA_A receptor data and on spike activity in an hiPSC-derived neuronal co-culture reported in this study are

Metric parameter	Description			
Mean spike rate (MSR)	Total number of spikes divided by recording time (Hz)			
ISI coefficient of variation	Standard deviation ISI (time between spikes) divided by the mean ISI. Measure for spike regularity: 0 indicates perfect spike distribution, >1 signals bursting			
Mean burst rate (MBR)	Total number of bursts divided by recording time (Hz)			
Burst duration	Average time from the first spike in a burst till the last spike (s)			
Number of spikes per burst	Average number of spikes occurring in a burst			
Mean inter-spike interval (ISI) within burst	Mean inter-spike interval within a burst (s)			
Inter-burst interval (IBI)	Time between the last spike of a burst and the first spike of a subsequent burst (s)			
IBI coefficient of variation	Standard deviation of IBI divided by the mean IBI. Measure for burst regularity			
Burst percentage	Percentage of total number of spikes occurring in a burst			
Mean network burst rate (MNBR)	Total number of network bursts divided by recording time (Hz)			
Network burst duration	Average time from the first spike till the last spike in a network burst (s)			
Number of spikes per network burst	Average number of spikes occurring in a network burst			
Mean ISI within network burst	Average of the mean ISIs within a network burst (s)			
Number of spikes per network burst per channel	Average number of spikes across the network bursts in the network burst divided by the number of electrodes participating in the network burst			
Network burst percentage	Percentage of total spikes occurring in a network burst			
Network IBI coefficient of variation	Standard deviation of network IBI divided by the mean network IBI. Measure of network burst rhythmicity: value is small when bursts occur at regular interval ar increases when bursts occur more sporadic			
Network normalized duration IQR	Interquartile range of network bursts durations. Measure for network burst duration regularity: larger values indicate wide variation in duration.			
Area under normalized cross-correlation	Area under inter-electrode cross-correlation normalized to the auto-correlations. The higher the value, the greater the synchronicity of the network			
Full width at half height (FWHH) of normalized cross-correlation	Width at half height of the normalized cross-correlogram. Measure for network synchrony: the higher the value, the less synchronized the network is			

Table 3. Description of different metric parameters. Adapted from⁵⁶.

within or below the range found in human plasma levels. It should be noted though that human exposure to PFAS is low dose, but continuous. Our *in vitro* study should therefore be considered mainly as hazard characterisation, highlighting the potential target(s) of PFOS and PFOA. Our findings, combined with the fact that PFAS can cross the blood-brain barrier, indicate that these compounds have a clear potential to cause neurotoxic effects in the occupationally exposed populations as well as in the general population.

Material and Methods

Animals. All experiments were conducted in accordance with Dutch law and approved by the Ethical Committee for Animal Experiments of Utrecht University. Animals were treated humanely and all efforts were made to alleviate suffering. *Xenopus laevis* oocytes were purchased from EcoCyte Bioscience (Castrop-Rauxel, Germany). Primary cultures of rat cortical neurons were prepared from postnatal day (PND) 0–1 Wistar rat pups (Envigo, Horst, The Netherlands).

Chemicals. MgSO₄, NaHCO₃ and Ca(NO₃) were purchased from Merck (Darmstadt, Germany). Neurobasal-A medium, Dulbecco's Modified Eagle Medium (DMEM)-F12, penicillin – streptomycin (5000 U/mL – 5000 µg/mL for rat primary cortical culture media and 10.000 U/mL – 10.000 µg/mL for hiPSC medium), B27 plus supplement, N2 supplement and L-glutamine were obtained from Life Technologies (Bleiswijk, The Netherlands). iCell Neural Supplement B and Nervous System Supplement were provided by Cellular Dynamics International (Madison, WI, USA). BrainPhys neuronal medium was obtained from StemCell Technologies (Cologne, Germany). SynFire seeding basal medium, short-term basal medium, long-term basal medium and accessory supplements were provided by NeuCyte (Sunnyvale, CA, USA). Laminin (L2020), 50% polyeth-yleneimine (PEI) solution, sodium borate, boric acid and all other chemicals (unless stated otherwise) were obtained from Sigma-Aldrich (Zwijndrecht, The Netherlands).

Stock solutions of PFOS (>98% purity, Fluka, Zwijndrecht, The Netherlands) and PFOA (>96% purity, Acros Organics, Geel, Belgium) were made in purity-checked dimethyl sulfoxide (DMSO) and stored at 4 °C. Final concentration of DMSO in PFOS- or PFOA-containing solutions was always kept below 0.1% (vol/vol). DMSO at concentrations up to 0.5% (vol/vol) had no effect on GABA_A receptor-mediated currents.

 $\alpha_1\beta_2\gamma_{2L}$ **GABA**_A receptor expression in Xenopus laevis oocytes. All procedures have been described previously^{43,48}. Briefly, complementary DNA (cDNA) coding for the human α_1 , β_2 (Origene, Rockville, USA) and human γ_{2L} subunits (kindly provided by Dr Paul J. Whiting, Merck Sharp & Dohme Research Laboratories, Neuroscience Research Center, Harlow, Essex, U.K.) was dissolved in distilled water at a 1:1:1 molar ratio and injected (23 nL/oocyte, ~1 ng of each subunit) into the nuclei of stage V or VI oocytes using a Nanoject Automatic Oocyte Injector (Drummond, Broomall, PA). Following injection, oocytes were incubated at 21 °C in modified

Barth's solution containing (in mM) 88 NaCl, 1 KCl, 2.4 NaHCO₃, 0.3 Ca(NO₃)₂, 0.41 CaCl₂, 0.82 MgSO₄, 15 HEPES, and 10μ g/mL neomycin (pH 7.6 with NaOH).

Electrophysiological recordings of GABA_A **receptor currents.** Electrophysiological recordings were performed on oocytes following 4–6 days of incubation to ensure sufficient translation of injected cDNA and functional expression of $\alpha_1\beta_2\gamma_{2L}$ GABA_A receptors in the membrane. Ion currents associated with GABA_A receptor activity were measured with the two-electrode voltage-clamp technique using a Gene Clamp 500B amplifier (Axon Instruments, Union City, CA) with high-voltage output stage as described previously ^{43,48}. Recording microelectrodes (0.1–1 MΩ) were filled with KCl (3 M). Oocytes, placed in a custom-built Teflon oocyte recording chamber, were voltage-clamped at –60 mV and continuously superfused (~30 mL/min) with saline solution containing (in mM) 1 CaCl₂, 10 HEPES (pH 7.2 with NaOH), 2.5 KCl, 115 NaCl. Membrane currents were low-pass filtered (8-pole Bessel; 3 dB at 0.3 kHz), digitized (12 bits; 1024 samples per record), and recorded on disk using WinWCP (Strathclyde Electrophysiological Software, Glasgow, Scotland) for further analysis.

Aliquots of freshly thawed GABA stock solution in distilled water and PFOS or PFOA in DMSO were added to the saline immediately before the experiments. GABA-responsive oocytes were exposed to saline, saline containing GABA (34μ M or 1 mM), saline containing a PFAS or saline containing GABA in combination with a PFAS by switching the perfusate using a servomotor-operated valve. To determine whether oocytes were GABA responsive, oocytes were super-fused for 40 s with a saline solution containing GABA. If oocytes were responsive, they were exposed for 10 s to saline containing GABA, followed directly by a 20 s exposure of saline containing GABA and PFAS and finally to 10 s of saline containing GABA. Concentrations of 0.01–100 μ M PFOS or PFOA were chosen to overlap with earlier *in vitro* studies and extended to include human serum and plasma concentrations. A washout period of 4 min between each application was introduced, allowing receptors to recover from desensitization. Sham-injected oocytes did not show any ion current upon superfusion with GABA, PFOS, or PFOA at concentrations tested, alone or in co-application with GABA (data not shown). To minimize adsorption of PFOS or PFOA to the perfusion system, glass reservoirs and Teflon tubes (polytetrafluoroethylene, Rubber, Hilversum, The Netherlands) were used.

Cell culture. Primary rat cortical cultures and hiPSC-derived neuronal co-cultures were kept at 37 °C in a humidified 5% CO_2 incubator. All cell culture surface materials were pre-coated with 0.1% PEI solution diluted in borate buffer (24 mM sodium borate/50 mM boric acid in Milli-Q adjusted to pH 8.4).

Primary rat cortical cells were isolated from PND0–1 Wistar rat pups as described previously^{50,64} with minor modifications. Briefly, PND0–1 pups were decapitated and cortices were rapidly dissected on ice and kept in serum free dissection medium (Neurobasal-A supplemented with 25 g/L sucrose, 450 μ M L-glutamine, 30 μ M glutamate, 1% penicillin/streptomycin and 2% B27 plus supplement, pH 7.4) during the entire procedure. Cortices were dissociated to a single-cell suspension by mincing with scissors, trituration and filtering through a 100 μ m mesh (EASY strainer, Greiner). The cell suspension was diluted to a 2 × 10⁶ cells/mL solution. Droplets of 50 μ L were placed on the electrode fields in wells of pre-coated 48-well MEA plates (Axion BioSystems Inc., Atlanta, GA, USA). Cells were left to adhere for ~2 hr before adding 450 μ L serum free dissection medium. At DIV4, 90% of the serum free dissection medium was replaced with serum free culture medium (Neurobasal-A supplemented with 25 g/L sucrose, 450 μ M L-glutamine, 1% penicillin/streptomycin and 2% B27 plus supplement, PH 7.4). Rat primary cortical neurons were exposed at DIV9–11.

iCell Glutaneurons (Lot# 103288, containing ~70% glutamatergic and ~30% GABAergic neurons); Cellular Dynamics International, Madison, WI, USA) and iCell Astrocytes (Lot# 103956; Cellular Dynamics International, Madison, WI, USA) were thawed and cultured according to manufacturer's protocol. In short, each cell type was thawed separately in supplemented BrainPhys medium (BrainPhys medium supplemented with 2% iCell Neural Supplement B, 1% Nervous System Supplement, 1% N2, 1% penicillin – streptomycin and 0.1% laminin). Following centrifugation, the cell pellet was diluted in dotting medium (supplemented BrainPhys medium with 10% laminin) till 15 × 10³ cells/µL for iCell Glutaneurons and 6.7 × 10³ cells/µL for iCell Astrocytes. Before plating, iCell Glutaneurons and iCell Astrocytes were premixed into a co-culture containing 120×10^3 iCell Glutaneurons and 20×10^3 iCell Astrocytes) over the electrode field of pre-coated 48-well MEA plates. After plating, cells were allowed to adhere for ~1 hr following which $300 \,\mu$ L of room temperature (RT) supplemented BrainPhys medium took place at DIV1, 2, 4, 6, 8, 10, 12 and 14. The iCell Glutaneurons – iCell Astrocytes cultures were exposed at DIV14.

SynFire glutamatergic neurons (Lot# 118B and 000124), SynFire GABAergic neurons (Lot# 118B and 000124) and SynFire astrocytes (Lot# 13029; all from NeuCyte Sunnyvalle, CA, USA) were thawed and cultured according to manufacturer's protocol. Briefly, each cell type was thawed separately in DMEM-F12. Each cell pellet was dissolved in complete seeding medium (containing the seeding supplement) till a density of 10×10^3 cells/µL following which a mixture was made containing 52% glutamatergic neurons, 22% GABAergic neurons and 26% astrocytes. This mixture was plated in 50µL droplets (270×10^3 cells per droplet) over the electrode field of pre-coated MEA plates. Cells were placed in the incubator and left overnight to adhere. The next day, 250µL/ well of complete short-term maintenance medium (containing the supplement) was added. At DIV3 and 5, 50% medium changes with short-term maintenance medium (containing supplements A and B) through 50% medium changes at DIV7, 10, 13, 16, 19, 22, 25 and 25. SynFire iNS co-cultures were exposed at DIV28.

MEA measurements. Each well of a 48-well MEA plate contains 16 nanotextured gold micro-electrodes (~40–50 µm diameter; 350 µm spacing) with 4 integrated ground electrodes yielding a total of 768 channels

for simultaneous recording (for review see⁷⁴). Spontaneous electrical activity was recorded as described previously^{52,64}. Briefly, signals were recorded at the day of experiments (DIV9–11 for rat primary cortical cultures, DIV14 for iCell Glutaneurons – iCell Astrocytes co-cultures and DIV28 for SynFire iNS co-cultures) using a Maestro 768-channel amplifier with integrated heating system and temperature controller and a data acquisition interface (Axion BioSystems Inc., Atlanta, GA, USA). Data acquisition was managed with Axion's Integrated Studio (AxIS 2.4.2.13) and recorded as.RAW files. All channels were sampled simultaneously with a gain of 1200x and a sampling frequency of 12.5 kHz/channel with a 200–5000 Hz band-pass filter. Prior to the recording, MEA plates were allowed to equilibrate for 5–10 min in the Maestro.

In order to determine effect of reference compounds on spontaneous neuronal activity (spiking and (network) bursting behaviour) on the three different cell cultures, a 30 min baseline recording was made. Following this baseline recording, wells were exposed ($10 \times$ dilution for primary rat cortical cultures and SynFire iNS co-culture and $30 \times$ dilution for iCell Glutaneurons – iCell Astrocytes co-culture) to PFOS, PFOA or the solvent control and directly a 30 min exposure recording was made. In order to overlap with earlier *in vitro* experiments and to fall in the range of reported human serum and plasma levels of PFOS and PFOA, concentrations of $0.1-100 \,\mu$ M were tested. Stock solutions of PFOS and PFOA in DMSO were diluted in culture medium to obtain desired concentrations. In all experiments, the solvent concentration never exceeded 0.1% v/v. In order to prevent receptor (de) sensitization each well was exposed to a single concentration. For each experimental condition, MEA plates from at least two different plating rounds were used.

Data analysis and statistics. Peak amplitudes of GABA-evoked ion currents were measured and normalized to the maximal amplitude (at 1 mM) of GABA-evoked control responses to adjust for differences in receptor expression levels among oocytes and for small variations in response amplitudes over time as described previously ^{43,48}. Normalized ion currents were plotted against GABA concentration in each experiment. GABA concentration-effect curves were fitted to the data obtained in separate experiments using Prism (Graphpad Software, La Jolla, CA, USA). The percentage of PFOS- or PFOA-induced inhibition of the GABA-evoked ion current was calculated from the quotient of the maximum amplitude of the GABA-PFOS/PFOA co-application response and the maximum amplitude of the control (GABA) response.

To determine (modulation of) spontaneous activity, RAW data files were re-recorded to obtain Alpha Map files. In this re-recording, spikes were detected with the AxIS spike detector (Adaptive threshold crossing, Ada BandFIt v2) using a variable threshold spike detector set at $7 \times$ (rat primary cortical cultures) or $5.5 \times$ (hiPSC cultures) standard deviation (SD) of internal noise level (rms) on each electrode. Post/pre-spike duration was set to 3.6/2.4 ms respectively. For further data analysis, spike files were loaded in NeuralMetric Tool (version 2.2.4, Axion BioSystems) and only active electrodes (MSR \geq 6 spikes/min) in active wells (\geq 1 active electrode) were included. Since (network) bursting behaviour is crucial for *in vivo* neuronal communication⁷⁵, it was analysed using the Poisson Surprise method⁷⁶ with a minimal surprise of 10 and a minimum bursting frequency of 0.005 bursts/sec. An adaptive threshold algorithm was used for the extraction of network bursts.

The effects of PFOS and PFOA on spontaneous neuronal activity were determined by comparing the baseline activity with activity following exposure. A custom-made MS Excel macro was used to calculate treatment ratios (TR) per well for the different metric parameters (Table 3) by: (parameter_{exposure}/parameter_{baseline}) × 100%. Hereafter, TRs were normalized to appropriate vehicle control (DMSO). To prevent inclusion of exposure artefacts, effect analysis was performed 20–30 min post-exposure.

Oocytes or wells that showed effect two times SD above or below average were considered outliers and removed for further data analysis (1.1% for oocytes, 4.2% for rat primary cortical neurons, 4.7% for iCell Glutaneuron – iCell Astrocytes co-cultures and 3.6% for SynFire iNS co-cultures). For oocyte data, the concentration-dependent effects of PFOS and PFOA were determined by one-way ANOVA and a post hoc Bonferroni test. Best-fit concentration-response curves were plotted in PRISM (version 8.0.1) by nonlinear regression with a four-parameter variable slope with the bottom value constrained to 0 and top value to 100 to improve the fit. In the case of MEA data, concentration-dependent effects were determined by a one-way Welch ANOVA and a post hoc Dunnett test. *P*-values < 0.05 were considered statistically significant. All statistical analyses were performed in R version 3.6.0 (R core team 2019) with base R or by using the DescTools package (by Signorell *et al.*, 2019, version 0.99.28).

Data are presented as mean \pm standard error of the mean (SEM) from the number of oocytes or wells (*n*) indicated, derived from 1 to 2 independent batches (*N*) of oocytes, 2 plating rounds (*N*) for the rat primary cortical culture, 5 plating rounds (*N*) for the iCell Glutaneuron – iCell Astrocytes co-culture and 2 independent plating rounds (*N*) for the SynFire iNS co-culture unless stated otherwise. IC₅₀ values are reported with confidence intervals (CI). LOECs are defined as lowest statistically significant concentrations.

Data availability

The datasets generated during and/or analysed during the current study are available from the corresponding author on reasonable request.

Received: 30 October 2019; Accepted: 4 March 2020; Published online: 24 March 2020

References

- 1. Grandjean, P. & Clapp, R. Perfluorinated Alkyl Substances: Emerging Insights Into Health Risks. *New Solut.* 25, 147–63 (2015).
- Lindstrom, A. B., Strynar, M. J. & Libelo, E. L. Polyfluorinated Compounds: Past, Present, and Future. *Environ. Sci. Technol.* 45, 7954–7961 (2011).

- 3. Environmental Protection Agency Federal Facilities Restoration, U. & Office, R. Technical Fact Sheet Perfluorooctane Sulfonate (PFOS) and Perfluorooctanoic Acid (PFOA). (2017).
- 4. Li, Y. *et al.* Half-lives of PFOS, PFHxS and PFOA after end of exposure to contaminated drinking water. *Occup. Environ. Med.* **75**, 46–51 (2018).
- 5. Olsen, G. W. *et al.* Half-life of serum elimination of perfluorooctanesulfonate, perfluorohexanesulfonate, and perfluorooctanoate in retired fluorochemical production workers. *Environ. Health Perspect.* **115**, 1298–305 (2007).
- Butenhoff, J. L., Ehresman, D. J., Chang, S.-C., Parker, G. A. & Stump, D. G. Gestational and lactational exposure to potassium perfluorooctanesulfonate (K+PFOS) in rats: Developmental neurotoxicity. *Reprod. Toxicol.* 27, 319–330 (2009).
- 7. Fuentes, S., Colomina, M. T., Vicens, P., Franco-Pons, N. & Domingo, J. L. Concurrent Exposure to Perfluorooctane Sulfonate and Restraint Stress during Pregnancy in Mice: Effects on Postnatal Development and Behavior of the Offspring. *Toxicol. Sci.* **98**, 589–598 (2007).
- Goulding, D. R., White, S. S., McBride, S. J., Fenton, S. E. & Harry, G. J. Gestational exposure to perfluorooctanoic acid (PFOA): Alterations in motor related behaviors. *Neurotoxicology* 58, 110–119 (2017).
- Johansson, N., Fredriksson, A. & Eriksson, P. Neonatal exposure to perfluorooctane sulfonate (PFOS) and perfluorooctanoic acid (PFOA) causes neurobehavioural defects in adult mice. *Neurotoxicology* 29, 160–9 (2008).
- 10. Onishchenko, N. *et al.* Prenatal exposure to PFOS or PFOA alters motor function in mice in a sex-related manner. *Neurotox. Res.* **19**, 452–61 (2011).
- 11. Wang, Y., Liu, W., Zhang, Q., Zhao, H. & Quan, X. Effects of developmental perfluorooctane sulfonate exposure on spatial learning and memory ability of rats and mechanism associated with synaptic plasticity. *Food Chem. Toxicol.* **76**, 70–76 (2015).
- 12. Pinkas, A., Ślotkin, T. A., Brick-Turin, Y., Van der Zee, E. A. & Yanai, J. Neurobehavioral teratogenicity of perfluorinated alkyls in an avian model. *Neurotoxicol. Teratol.* **32**, 182–186 (2010).
- Spulber, S. et al. PFOS Induces Behavioral Alterations, Including Spontaneous Hyperactivity That Is Corrected by Dexamfetamine in Zebrafish Larvae. PLoS One 9, e94227 (2014).
- Ulhaq, M., Orn, S., Carlsson, G., Morrison, D. A. & Norrgren, L. Locomotor behavior in zebrafish (Danio rerio) larvae exposed to perfluoroalkyl acids. Aquat. Toxicol. 144–145, 332–40 (2013).
- Liew, Z. et al. Prenatal Exposure to Perfluoroalkyl Substances and the Risk of Congenital Cerebral Palsy in Children. Am. J. Epidemiol. 180, 574–581 (2014).
- Høyer, B. B. et al. Pregnancy serum concentrations of perfluorinated alkyl substances and offspring behaviour and motor development at age 5-9 years - a prospective study. Environ. Heal. 14, 2 (2015).
- 17. Harris, M. H. *et al.* Prenatal and childhood exposure to per- and polyfluoroalkyl substances (PFASs) and child cognition. *Environ. Int.* **115**, 358–369 (2018).
- Forns, J. et al. Perfluoroalkyl substances measured in breast milk and child neuropsychological development in a Norwegian birth cohort study. Environ. Int. 83, 176–182 (2015).
- Liew, Z. et al. Prenatal Exposure to Perfluoroalkyl Substances and IQ Scores at Age 5; a Study in the Danish National Birth Cohort. Environ. Health Perspect. 126, 067004 (2018).
- Liew, Z. et al. Attention deficit/hyperactivity disorder and childhood autism in association with prenatal exposure to perfluoroalkyl substances: a nested case-control study in the Danish National Birth Cohort. Environ. Health Perspect. 123, 367–73 (2015).
- Strøm, M. et al. Persistent organic pollutants measured in maternal serum and offspring neurodevelopmental outcomes A
 prospective study with long-term follow-up. Environ. Int. 68, 41–48 (2014).
- Goudarzi, H. et al. Prenatal exposure to perfluorinated chemicals and neurodevelopment in early infancy: The Hokkaido Study. Sci. Total Environ. 541, 1002–1010 (2016).
- 23. Quaak, I. et al. Prenatal exposure to perfluoroalkyl substances and Behavioral development in children. Int. J. Environ. Res. Public Health 13, (2016).
- 24. Ghassabian, A. *et al.* Concentrations of perfluoroalkyl substances and bisphenol A in newborn dried blood spots and the association with child behavior. *Environ. Pollut.* **243**, 1629–1636 (2018).
- Lenters, V. et al. Early-life exposure to persistent organic pollutants (OCPs, PBDEs, PCBs, PFASs) and attention-deficit/hyperactivity disorder: A multi-pollutant analysis of a Norwegian birth cohort. Environ. Int. 125, 33–42 (2019).
- Berntsen, H. F. et al. Time-dependent effects of perfluorinated compounds on viability in cerebellar granule neurons: Dependence on carbon chain length and functional group attached. *Neurotoxicology* 63, 70–83 (2017).
- Chen, N., Li, J., Li, D., Yang, Y. & He, D. Chronic exposure to perfluorooctane sulfonate induces behavior defects and neurotoxicity through oxidative damages, *in vivo* and *in vitro*. *PLoS One* 9, e113453 (2014).
- Reistad, T., Fonnum, F. & Mariussen, E. Perfluoroalkylated compounds induce cell death and formation of reactive oxygen species in cultured cerebellar granule cells. *Toxicol. Lett.* 218, 56–60 (2013).
- 29. Zhang, L. et al. Perfluorooctane Sulfonate Induces Apoptosis in N9 Microglial Cell Line. Int. J. Toxicol. 30, 207-215 (2011).
- Slotkin, T. A., MacKillop, E. A., Melnick, R. L., Thayer, K. A. & Seidler, F. J. Developmental Neurotoxicity of Perfluorinated Chemicals Modeled in Vitro. Environ. Health Perspect. 116, 716–722 (2008).
- Lee, H.-G., Lee, Y. J. & Yang, J.-H. Perfluorooctane sulfonate induces apoptosis of cerebellar granule cells via a ROS-dependent protein kinase C signaling pathway. *Neurotoxicology* 33, 314–320 (2012).
- 32. Mashayekhi, V. et al. Mechanistic approach for the toxic effects of perfluorooctanoic acid on isolated rat liver and brain mitochondria. Hum. Exp. Toxicol. 34, 985–96 (2015).
- Sun, P. et al. Nrf2 Signaling Elicits a Neuroprotective Role Against PFOS-mediated Oxidative Damage and Apoptosis. Neurochem. Res. 43, 2446–2459 (2018).
- Wang, C. et al. Reactive oxygen species mediate nitric oxide production through ERK/JNK MAPK signaling in HAPI microglia after PFOS exposure. Toxicol. Appl. Pharmacol. 288, 143–151 (2015).
- Liu, X., Jin, Y., Liu, W., Wang, F. & Hao, S. Possible mechanism of perfluorooctane sulfonate and perfluorooctanoate on the release of calcium ion from calcium stores in primary cultures of rat hippocampal neurons. *Toxicol. Vitr.* 25, 1294–1301 (2011).
- 36. Yang, J. *et al.* Perfluorooctane sulfonate mediates microglial activation and secretion of TNF-α through Ca2+-dependent PKC-NFκB signaling. *Int. Immunopharmacol.* **28**, 52–60 (2015).
- Liao, C.-Y., Cui, L., Zhou, Q.-F., Duan, S.-M. & Jiang, G.-B. Effects of perfluorooctane sulfonate on ion channels and glutamateactivated current in cultured rat hippocampal neurons. *Environ. Toxicol. Pharmacol.* 27, 338–44 (2009).
- Liao, C. et al. Changes in synaptic transmission, calcium current, and neurite growth by perfluorinated compounds are dependent on the chain length and functional group. Environ. Sci. Technol. 43, 2099–104 (2009).
- Liao, C.-Y., Li, X.-Y., Wu, B., Duan, S. & Jiang, G.-B. Acute enhancement of synaptic transmission and chronic inhibition of synaptogenesis induced by perfluorooctane sulfonate through mediation of voltage-dependent calcium channel. *Environ. Sci. Technol.* 42, 5335–41 (2008).
- 40. Zhang, Q. *et al.* Developmental perfluorooctane sulfonate exposure inhibits long-term potentiation by affecting AMPA receptor trafficking. *Toxicology* **412**, 55–62 (2019).
- Antunes Fernandes, E. C., Hendriks, H. S., van Kleef, R. G. D. M., van den Berg, M. & Westerink, R. H. S. Potentiation of the Human GABA A Receptor As a Novel Mode of Action of Lower-Chlorinated Non-Dioxin-Like PCBs[†]. *Environ. Sci. Technol.* 44, 2864–2869 (2010).

- 42. Antunes Fernandes, E. C. *et al.* Activation and Potentiation of Human GABAA Receptors by Non-Dioxin–Like PCBs Depends on Chlorination Pattern. *Toxicol. Sci.* **118**, 183–190 (2010).
- Hendriks, H. S., Antunes Fernandes, E. C., Bergman, Å., van den Berg, M. & Westerink, R. H. S. PCB-47, PBDE-47, and 6-OH-PBDE-47 Differentially Modulate Human GABAA and α4β2 Nicotinic Acetylcholine Receptors. *Toxicol. Sci.* 118, 635–642 (2010).
- Hendriks, H. S., van Kleef, R. G. D. M., van den Berg, M. & Westerink, R. H. S. Multiple Novel Modes of Action Involved in the *In Vitro* Neurotoxic Effects of Tetrabromobisphenol-A. *Toxicol. Sci.* 128, 235–246 (2012).
- 45. Represa, A. & Ben-Ari, Y. Trophic actions of GABA on neuronal development. Trends Neurosci. 28, 278-283 (2005).
- Fujii, S., Jia, Y., Yang, A. & Sumikawa, K. Nicotine reverses GABAergic inhibition of long-term potentiation induction in the hippocampal CA1 region. Brain Res. 863, 259–65 (2000).
- Olsen, R. W. & Sieghart, W. GABAA receptors: Subtypes provide diversity of function and pharmacology. *Neuropharmacology* 56, 141–148 (2009).
- Hondebrink, L. et al. Structure-dependent inhibition of the human α1β2γ2 GABAA receptor by piperazine derivatives: A novel mode of action. Neurotoxicology 51, 1–9 (2015).
- 49. Robinette, B. L., Harrill, Ja, Mundy, W. R. & Shafer, T. J. *In Vitro* Assessment of Developmental Neurotoxicity: Use of Microelectrode Arrays to Measure Functional Changes in Neuronal Network Ontogeny1. *Front. Neuroeng.* **4**, 1 (2011).
- Dingemans, M. M. L. et al. Chronic 14-day exposure to insecticides or methylmercury modulates neuronal activity in primary rat cortical cultures. Neurotoxicology 57, 194–202 (2016).
- McConnell, E. R., McClain, M. A., Ross, J., LeFew, W. R. & Shafer, T. J. Evaluation of multi-well microelectrode arrays for neurotoxicity screening using a chemical training set. *Neurotoxicology* 33, 1048–1057 (2012).
- 52. Nicolas, J. et al. Detection of marine neurotoxins in food safety testing using a multielectrode array. Mol. Nutr. Food Res. 58, 2369–2378 (2014).
- Vassallo, A. et al. A multi-laboratory evaluation of microelectrode array-based measurements of neural network activity for acute neurotoxicity testing. Neurotoxicology 60, 280–292 (2017).
- Little, D., Ketteler, R., Gissen, P. & Devine, M. J. Using stem cell-derived neurons in drug screening for neurological diseases. *Neurobiol. Aging* 78, 130–141 (2019).
- Tukker, A. M., Wijnolts, F. M. J., de Groot, A. & Westerink, R. H. S. Human iPSC-derived neuronal models for *in vitro* neurotoxicity assessment. *Neurotoxicology* 67, 215–225 (2018).
- Tukker, A. M., Van Kleef, R. G. D. M., Wijnolts, F. M. J., De Groot, A. & Westerink, R. H. S. Towards animal-free neurotoxicity screening: Applicability of hiPSC-derived neuronal models for *in vitro* seizure liability assessment. *ALTEX* 37, 121–135 (2020).
- Mariussen, E. Neurotoxic effects of perfluoroalkylated compounds: mechanisms of action and environmental relevance. Arch. Toxicol. 86, 1349–1367 (2012).
- D'Hulst, C., Atack, J. R. & Kooy, R. F. The complexity of the GABAA receptor shapes unique pharmacological profiles. *Drug Discov. Today* 14, 866–875 (2009).
- Berntsen, H. F. et al. PFOS-induced excitotoxicity is dependent on Ca2+ influx via NMDA receptors in rat cerebellar granule neurons. Toxicol. Appl. Pharmacol. 357, 19–32 (2018).
- Ishida, K., Tsuyama, Y., Sanoh, S., Ohta, S. & Kotake, Y. Perfluorooctane sulfonate induces neuronal vulnerability by decreasing GluR2 expression. Arch. Toxicol. 91, 885–895 (2017).
- 61. Neucyte Inc. NeuCyte|Data. Available at: https://www.neucyte.com/data. (Accessed: 11th September 2019).
- International, C. D. iCell GlutaNeurons|Cellular Dynamics. Available at: https://fujifilmcdi.com/products-services/icell-products/ glutaneurons/. (Accessed: 3rd January 2020)
- Görtz, P., Fleischer, W., Rosenbaum, C., Otto, F. & Siebler, M. Neuronal network properties of human teratocarcinoma cell linederived neurons. *Brain Res.* 1018, 18–25 (2004).
- 64. Tukker, A. M. *et al.* Is the time right for *in vitro* neurotoxicity testing using human iPSC-derived neurons? *ALTEX* **33**, 261–271 (2016).
- 65. Tang, X. *et al*. Astroglial cells regulate the developmental timeline of human neurons differentiated from induced pluripotent stem cells. *Stem Cell Res.* **11**, 743–757 (2013).
- Ishii, M. N., Yamamoto, K., Shoji, M., Asami, A. & Kawamata, Y. Human induced pluripotent stem cell (hiPSC)-derived neurons respond to convulsant drugs when co-cultured with hiPSC-derived astrocytes. *Toxicology* 389, 130–138 (2017).
- 67. Wu, X. *et al.* Astrocytes are protective against chlorpyrifos developmental neurotoxicity in human pluripotent stem cell-derived astrocyte-neuron cocultures. *Toxicol. Sci.* **157**, 410–420 (2017).
- Li, Z. et al. Evaluation of PFOS-mediated neurotoxicity in rat primary neurons and astrocytes cultured separately or in co-culture. Toxicol. Vitr. 38, 77–90 (2017).
- Harada, K. H., Ishii, T. M., Takatsuka, K., Koizumi, A. & Ohmori, H. Effects of perfluorooctane sulfonate on action potentials and currents in cultured rat cerebellar Purkinje cells. *Biochem. Biophys. Res. Commun.* 351, 240–245 (2006).
- Ehresman, D. J., Froehlich, J. W., Olsen, G. W., Chang, S.-C. & Butenhoff, J. L. Comparison of human whole blood, plasma, and serum matrices for the determination of perfluorooctanesulfonate (PFOS), perfluorooctanoate (PFOA), and other fluorochemicals. *Environ. Res.* 103, 176–184 (2007).
- Fromme, H., Tittlemier, S. A., Völkel, W., Wilhelm, M. & Twardella, D. Perfluorinated compounds Exposure assessment for the general population in western countries. Int. J. Hyg. Environ. Health 212, 239–270 (2009).
- Wang, J. et al. Penetration of PFASs Across the Blood Cerebrospinal Fluid Barrier and Its Determinants in Humans. Environ. Sci. Technol. 52, 13553–13561 (2018).
- 73. Wu, M. et al. Analysis of perfluorinated compounds in human serum from the general population in Shanghai by liquid chromatography-tandem mass spectrometry (LC-MS/MS). Chemosphere 168, 100-105 (2017).
- Johnstone, A. F. M. et al. Microelectrode arrays: A physiologically based neurotoxicity testing platform for the 21st century. NeuroToxicology 31, 331–350 (2010).
- Izhikevich, E. M., Desai, N. S., Walcott, E. C. & Hoppensteadt, F. C. Bursts as a unit of neural information: selective communication via resonance. *Trends Neurosci.* 26, 161–167 (2003).
- Legéndy, C. R. & Salcman, M. Bursts and recurrences of bursts in the spike trains of spontaneously active striate cortex neurons. J. Neurophysiol. 53, 926–939 (1985).

Acknowledgements

This work was funded by a grant from the National Centre for the Replacement, Refinement and Reduction of Animals in Research (NC3Rs; project number 50308–372160). Members of the Neurotoxicology Research Group are gratefully acknowledged for helpful discussions. We thank Fiona Wijnolts for her assistance with cell culture work and Barbara Vreede (University library, Utrecht University) for her help with R scripts.

Author contributions

All authors contributed to the study conception and design. Material preparation, data collection and analysis were performed by A.T., L.B. and G.K. The first draft of the manuscript was written by A.T. and H.H. R.W. and J.L. were involved with study design and helpful discussion concerning the results and manuscript. All authors commented on previous versions of the manuscript. All authors read and approved the final manuscript.

Competing interests

The authors declare no competing interests.

Additional information

Supplementary information is available for this paper at https://doi.org/10.1038/s41598-020-62152-2.

Correspondence and requests for materials should be addressed to R.H.S.W.

Reprints and permissions information is available at www.nature.com/reprints.

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

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 license, and indicate if changes were made. The images or other third party material in this article are included in the article's Creative Commons license, unless indicated otherwise in a credit line to the material. If material is not included in the article's Creative Commons license 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 license, visit http://creativecommons.org/licenses/by/4.0/.

© The Author(s) 2020