

SCIENCE COPYWRITER

Why Axion BioSystems?

At Axion, we empower our scientists and engineers to develop the tools that lead to new discoveries and better clinical outcomes. Our products are actively used in ground-breaking research in the fields of neurodegenerative disease, cancer immunotherapy, and cardiac safety. Located in Atlanta, GA, you'll join a team that is skilled, passionate, and supportive. If you want to be a part of a company that is on the leading edge of scientific research, Axion may be the place for you.

Job Description Summary:

We are looking for a Science Copywriter to join our marketing team at Axion BioSystems. The Science Copywriter is responsible for producing clear, concise and engaging content for web, social media, online publications and print materials to support marketing projects for Axion BioSystems. The right candidate must have the ability to manage multiple projects and deadlines and prioritize tasks in a fast-paced environment. They will be capable of translating scientific research into engaging pieces that can be easily read by Axion's audience. They can adapt their writing style to ensure all messaging and copy is within brand voice and editorial guidelines and are capable of writing long- or short-form copy.

Required Skills:

- Bachelor's degree in Communication or related field or an advanced degree in biological sciences with demonstrated skill in writing.
- 3-5 years in a professional copywrite role
- Skilled at translating scientific material to a consumer audience
- Aptitude for managing multiple projects, changing deadlines, and working in teams.
- Excellent verbal, written, and interpersonal communication skills.
- Ability to work both independently of and cooperatively with team members.
- Writing should be stellar in grammar, style and tone. Short and long-form copy should be versatile, as well as effective for headlines, taglines, and calls to action.
- A thorough knowledge of contemporary editorial and web practices.
- Fluent in written and spoken English
- U.S Citizenship or permanent residency required

Preferred Skills:

The following are not required, but will enhance an applicant's candidacy.

- Experience in biotech or biology-related field strongly preferred

Please provide examples of previous work.

CONTACT INFORMATION:

If you are interested in this position, please send your resume and cover letter to careers@axionbio.com

