

## SOCIAL MEDIA SPECIALIST

### 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 creative Social Media Specialist to join our Marketing team at Axion BioSystems. As the Social Media Specialist, you will work with our digital marketing team to manage Axion's social media presence. You will develop, schedule and maintain Axion's corporate social media channels and engage with our online scientific community. Responsibilities will include: developing content that resonates with our scientific audience; implement and monitor campaigns that align with Axion initiatives; collect, report and manage social media metrics from both organic and paid analytics. The ideal candidate will have previous experience managing social media marketing campaigns and share our passion for communicating the latest in scientific research.

### Required Skills:

- Bachelor of Science in Biology, Biomedical Engineering, or similar field
- 1-3 years of work experience in content marketing or communications
- Comprehend scientific writing at a high level of science and encapsulate the findings
- Attention to detail with excellent communication skills
- Fluent in written and spoken English
- Aptitude for managing multiple projects, changing deadlines, and working in teams
- U.S Citizenship or permanent residency required

### Preferred Skills:

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

- Advanced experience with online content searches and current knowledge of internet trends in scientific context
- Demonstrate success in social media marketing, including social media metrics
- Proficiency with multiple social media platforms such as LinkedIn, YouTube, and Twitter.

### CONTACT INFORMATION:

If you are interested in this position, please send your resume and cover letter to [careers@axionbio.com](mailto:careers@axionbio.com)

