

## MARKETING GRAPHIC DESIGNER

### 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 Marketing Graphic Designer to join our Marketing team at Axion BioSystems. As the Marketing Graphic Designer, you will support creative strategy and design/visual standards to develop superior creative work across all department deliverables. Responsibilities will include: design and manage Axion brand assets and customer-facing collateral, ensure consistent visual standards and content proofing, produce and edit promotional videos, coordinate with our digital team to develop graphics and website assets, collaborate with internal and external scientists for the production of marketing material. The ideal candidate will have good communication, creativity, an attention to detail, and share our passion for communicating the latest in scientific research.

### Required Skills:

- Associate's degree in design, marketing, or biological sciences. Bachelor's preferred
- 1-3 years of marketing or graphic design experience
- Aptitude for managing multiple projects, changing deadlines, and working in teams.
- Intermediate proficiency in Adobe Illustrator, Adobe Photoshop, and Adobe Premiere
- Experience with Adobe InDesign, Adobe After Effects
- Must display, by portfolio, superior design and art direction skills
- Must be a creative, conceptual thinker
- Attention to detail with excellent communication skills
- 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 with photography, videography, and video editing
- Experience in biotech or biology-related field strongly preferred

**Please provide a link to your portfolio or examples of previous work.**

### CONTACT INFORMATION:

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

