



1914 Palomar Oaks Way
Suite 150
Carlsbad CA 92008

WEB CONTENT COORDINATOR

Active Motif, Inc. is a privately held biotechnology company focused on supplying innovative kits and reagents for epigenetics and nuclear function research. Active Motif's Marketing group is currently recruiting highly motivated and talented candidates for a Web Content Coordinator position to join our team.

This role will be responsible for working with Product Managers to design content for web and email campaigns. A successful candidate will work closely with other members of the IT and marketing teams and communicate progress and challenges to management. This is an excellent opportunity for candidates seeking to apply self-taught/freelance skills in a professional environment or for skilled candidates looking to take a step forward in their career.

Job Responsibilities:

- Gather and evaluate user requirements and develop user stories in collaboration with product managers and developers.
- Illustrate design ideas using storyboards, process flows, and sitemaps.
- Develop low and high fidelity UI mockups and prototypes that clearly illustrate function and feel.
- Manage multiple projects effectively and efficiently while maintaining priorities, deadlines, and deliverables.
- Communicate issues and concerns to stakeholders and leadership frequently and clearly.
- Design HTML layouts for web & email content.
- Create banners, images, promotions, etc. for email marketing (and other marketing) campaigns on Active Motif and/or 3rd party vendor websites, using Adobe's Creative Suite (Photoshop, Illustrator, InDesign, etc.)

Required Qualifications:

- Outstanding HTML/CSS skills.
- Experience creating and populating web content in one or more commonly-used Content Management Systems using a WYSIWYG editor, as well as in HTML code.
- Relational database querying with SQL. Familiarity with a GUI database management tool (eg. HeidiSQL, MySQL Workbench) is a plus.
- Experience with Adobe Creative Cloud suite (Photoshop, Illustrator, InDesign, etc.).
- Thorough understanding of basic web technologies, frameworks and concepts (FTP, CMS, SEO, analytics, e-commerce, web design patterns, etc.) as they relate to content design and management.
- Strong technical troubleshooting skills.
- Understanding of PHP a plus.
- A background or experience in molecular biology would be useful.

In addition to a highly interactive and stimulating work environment, Active Motif also provides competitive salaries, stock options, and an excellent benefits package. Active Motif is an equal opportunity employer and promotes diversity throughout its workforce. Active Motif is located in Carlsbad, California, and has offices in Belgium, Japan and China.

If you are interested in a career with Active Motif, please submit your cover letter and resume outlining qualifications, experience and current salary details to hr@activemotif.com. We thank all applicants, however, only those selected for interviews will be contacted.