



1 Great Valley Parkway, Suite 2, Malvern, PA 19355, USA
1-610-722-0247 • www.reactionbiology.com

Title: Scientific Marketing Communications Manager

Job description

Reaction Biology is a private contract research organization supporting the drug discovery programs of over 1800 clients worldwide from pharma, biotech or academia through our operations in Malvern, PA and Freiburg, Germany.

We are looking for a scientific marketing communications manager to promote the Reaction Biology portfolio and brand via execution of marketing campaigns in a range of offline and online marketing channels. The candidate should bring marketing experience and a science background.

Job responsibilities

- Implement marketing communication plans by coordination of diverse marketing activities and execution of marketing campaigns
- Maximize marketing visibility and promotion of our brand values across all forms of media and advertisement outlets
- Coordinate relationship with publishers and online platforms
- Work closely with the marketing director to create marketing programs
- Generate print and online creatives in cooperation with a designer
- Coordinate printing and dissemination of print marketing material
- Manage social media presence, display advertisement, newsletters and email messaging
- Collect metrics and customer behavioral data in response to marketing activities
- Prepare competitive market analysis to identify strategic options to grow market share
- Plan and organize the logistics of our presence at conferences and trade shows including booth assembly and coordination of marketing activities

Position qualifications

- Required: Scientific education and at least 2 years of marketing-related experience in a biotechnology or pharmaceutical company
- Knowledge of common marketing communication practices
- Demonstrated initiative in implementation of marketing communication plans
- Excellent written communication skills ideally demonstrated interest in creative writing
- High level of organization and attention to detail
- Collaborative and proactive attitude
- Ability to work independently and as part of a team

Let's discover together.



- Flexibility to reprioritize quickly to meet changing priorities and deliver results on deadline
- Asset: Knowledge of the Adobe creative suite

Reporting to the Marketing Director

Employment type

- Full time
- Location: Remote or at our location in Malvern, PA, USA
- Citizenship: any

