



**Job Posting: Lab Assistant**

Echelon Biosciences is looking for a part-time lab assistant to join our team. Echelon is a Salt Lake City-based, world-recognized supplier of novel research and drug discovery products focused on cell signaling and diagnostic reagents. We research, develop, and manufacture reagents/assay products and provide excellent technical support for academic, biotech, and pharmaceutical customers.

**Position Requirements:**

- A high school diploma or equivalent
- Experience working in a scientific lab or biotechnology company a plus
- Ability to work both independently as well as part of a team
- Effective written and oral communication and organizational skills
- Part-time position (12 hours per week)

**Position Tasks and Responsibilities:**

- Weigh out of finished goods product for orders (includes weighing material, labeling vials, packaging orders)
- Inventory management (proper cataloging of vial containers, low/no quantity tracking, quality control tracking, backorder tracking/communication)
- Coordinate with other personnel regarding inventory needs and shipping of orders
- Help with tasks in lab and office as needed
- Miscellaneous projects throughout year (as needed) such as year-end inventory count

For more information, visit our website at [www.echelon-inc.com](http://www.echelon-inc.com), email your resume to [echelon@echelon-inc.com](mailto:echelon@echelon-inc.com) or mail to 675 Arapeen Drive Ste 302 Salt Lake City, UT 84108.

Echelon Biosciences, Inc. is an Equal Opportunity Employer.