
Request for Applications – Biotechnology Internship, Summer 2023

Echelon Biosciences is a biologic reagents company based in Salt Lake City, UT. We are seeking applications for our Internship Program from qualified and highly motivated university students that are interested in pursuing a career in biotechnology and related fields.

Our program's goal is to encourage young, aspiring scientists by providing them with a unique environment within the biotechnology sector for project-based training.

Successful applicants will be assigned a project mentor based on the applicants' stated goals and knowledge base. Successful applicants may be asked to assist with a variety of activities ranging from basic product quality procedures to product research and development.

General application guidelines are as follows:

- Applicants must be enrolled in a relevant degree program related to the **Biological Sciences, Bioengineering, or Chemistry**
- Applicants must be able to commit to a specified work load of at least half-time = 15 hrs/wk for 12 weeks
- Must have completed science course pre-requisites and be enrolled in upper-level courses; minimum two years of completed course work in total
- Applicants will also be expected to have good organizational and time management skills

Start dates for individual internships are also flexible based on the applicant's university schedule.

If interested, please send a resumé with relevant course work, grades, at least one reference, and any additional experience along with a cover letter to: echelon@echelon-inc.com, attn: Internship Program. Resumés and applications may also be submitted via [Handshake.com](https://www.handshake.com).

Anticipated start date for the summer term is **June 1**. We are currently recruiting for the following projects:

Characterization of Lipid Antibodies for Immunostaining

Project seeks to characterize a panel of antibodies for specific lipids under different chemical staining conditions.

Successful applicants will gain relevant experience with cell culture, immunocytochemistry, antibody handling, and protein purification.

Development of Lateral Flow/In Vitro Diagnostic for Alcohol Biomarker

Project is working toward transfer of existing product technology to a point-of-care based assay for a known alcohol biomarker.

Successful applicants will gain relevant experience with lipid preparation and extraction, handling synthetic and human samples, and clinical assay development.