



## RESEARCH TRAVEL GRANT APPLICATION

Enzo Research Travel Grant is available for scientists who are presenting their work featuring Enzo's products at research conferences and meetings. Enzo will provide up to \$1,000 to cover travel expenses.

---

**Name:** \_\_\_\_\_ **Degree/Title:** \_\_\_\_\_

**Email Address:** \_\_\_\_\_

**Phone Number:** \_\_\_\_\_

**Institution:** \_\_\_\_\_

**Principal Investigator:** \_\_\_\_\_

**Job Title:** \_\_\_\_\_

**Department:** \_\_\_\_\_

**Address:** \_\_\_\_\_

**City:** \_\_\_\_\_ **State:** \_\_\_\_\_ **Zip:** \_\_\_\_\_

**Country:** \_\_\_\_\_



**Conference:** \_\_\_\_\_

**Dates:** \_\_\_\_\_

**Location:** \_\_\_\_\_

**Project Title:** \_\_\_\_\_

**Product(s) Featured:**

**Abstract**



**ABOUT ENZO RESEARCH TRAVEL GRANT:** Enzo Biochem is offering grant awards for scientists to help defray travel-related expenses associated with attending conferences and meetings. These funds are meant to defray the cost of lodging and/or air/ground transportation expenses. Scientists presenting work featuring Enzo's products are eligible to apply for the travel grants. The amount of the award is up to \$1,000 maximum and must be applied to related expenses, including travel (air and ground transportation) and meeting registration.

**These awards will be processed as reimbursements to the award recipients after the meeting is over. Original receipts will be required in order to receive reimbursement. The maximum amount to be reimbursed for meeting- and/or travel-related expenses will be indicated in the award recipient's award letter.**

If you apply and are approved to receive a travel grant and later change your mind about presenting your work, the entire grant will be rescinded and cancelled completely, and you will not be reimbursed for any expenses you have already incurred, or eventually incur.

**By checking this box, you acknowledge that you read, understand, and fully accept the terms and conditions for this award. You, the applicant, represent that you understand the above terms and conditions as written in English or had the terms and conditions translated into the appropriate language in order to comprehend the terms and conditions for this award. To the extent you translated the aforementioned terms and conditions, you fully assume all risks associated with translating these terms and conditions.**

ELIGIBLE SCIENTIST WHO WISH TO BE CONSIDERED FOR ENZO TRAVEL GRANT MUST:

1. Follow abstract submission guidelines of the conference you are attending. When you receive your notification letter indicating that your research abstract will be programmed and included as a poster or oral presentation at the conference or meeting, you may complete the Enzo Travel Grant application process by submitting the confirmation letter ASAP, and no later than 14 calendar days before the first day of the conference. Failure to complete this process will void your Enzo Research Travel Grant application.
2. Complete and submit this application form including all supporting documentation.
  - a. A copy of your **letter of confirmation** of programmed oral/poster presentation (within 48 hours after notification of your confirmed programmed abstract; no later than 14 calendar days before the first day of the meeting.)
  - b. A copy of your current CV/resume.

All grant applicants must be approved by Enzo's Grant review committee. Abstract submitted and accepted as poster or platform presentations will be used as the basis for the awards. Enzo's Grant review committee will select the award recipients based on all supporting documentations from applicants who are eligible. Funding for the awards are contingent upon funding availability. Applications will be reviewed and approved on a rolling basis.

Applicants selected to receive an award will be notified by Enzo Grant program staff upon approval of their application.

SUBMIT APPLICATION TO: [grants@enzo.com](mailto:grants@enzo.com).



## ENZO RESEARCH TRAVEL GRANT APPLICATION CHECKLIST

- Complete travel grant application
- Copy of the confirmation/notification of abstract acceptance for meeting program
- Copy of CV/Resume

**Please be advised:** We do not accept piecemeal submissions. All incomplete or late application packets will be declined upon receipt.

## TERMS AND CONDITIONS

By submitting your entry, you agree to these terms for yourself, on behalf of the principal investigator (PI) and your Institution. “You” and “your” shall refer to you personally, your PI, or your Institution, as applicable. “We,” “Us,” or “Our” includes Enzo Biochem and/or its officers, employees, directors, agents, and affiliates. If you do not agree to these terms, do not participate in this offer.

1. Only one entry per individual. Once entry is submitted, you may not amend it. Applications must be submitted at least 14 calendar days prior to the first day of the conference to qualify.
2. All qualified entries will be judged by a team of Enzo scientists and grant review committee for appropriateness and completeness in response.
3. Successful grant applicants will receive up to \$1,000 to help defray travel expenses incurred to attend the conference where an oral and/or poster presentation is programmed.
4. Grant award winners agree to make their project datasets available for confidential internal review by Enzo within 120 days following notification of winning an award.
5. Grant winners agree to be available for reasonable promotional activities at Enzo’s request for a period of twelve months following announcement of a prize award. Promotional activities may include, but are not limited to:
  - Assisting Enzo in preparation of a case study or application note based on the project
  - Participating in Enzo’s webinars
  - Being featured as a researcher in an Enzo Spotlight/Website/Social Media story
  - Participating in virtual or in-person seminars
  - Image release
6. You agree to obtain all authorizations and permissions from your Institution and PI or others that are required to allow you to make this submission and to have your name and the name of your institution used for all purposes, including marketing, by Enzo or its affiliates without any compensation and with or without attribution, regardless of whether yours is judged to be a winning entry. You also agree personally to all the authorizations and permissions required by these terms and conditions.
7. You agree that the use of your submission by Enzo does not violate any rule, policy, copyright, trademark, patent, or procedures of your Institution or that of any other person. Participation in submission does not implead on intellectual property of the results.
8. You acknowledge and agree that any submission made hereunder is not being made in exchange for any promise of providing business to Enzo.



9. You agree that Enzo may contact you for purposes of marketing its products and services.
10. Enzo reserves the right to discontinue or modify this program at any time for any reason. In such event, Enzo will notify and inform all already-submitted entrants via e-mail. In the event of discontinuance or modification of the program, this will not result in Enzo having any liability to you, your employees, agents, or team members for such discontinuance or modification.
11. Enzo may refuse to consider a submission if Enzo determines in its sole discretion that awarding the prize to the entrant would likely result in a violation of an applicable law, rule, or policy, including, but not limited to, whether awarding a prize would be contrary to the law where you reside or work.
12. IN NO EVENT SHALL ENZO, ITS OFFICERS, DIRECTORS, EMPLOYEES, OR AGENTS, BE LIABLE FOR ANY DIRECT, INDIRECT, INCIDENTAL, SPECIAL, PUNITIVE, OR CONSEQUENTIAL DAMAGES WHATSOEVER ARISING UNDER OR AS A RESULT OF ENTRANCE IN OR PARTICIPATION IN THE OFFER, WHETHER BASED ON WARRANTY, CONTRACT, TORT, STRICT LIABILITY, OR ANY OTHER LEGAL THEORY, AND WHETHER OR NOT ENZO IS ADVISED OF THE POSSIBILITY OF SUCH DAMAGES. THE FOREGOING LIMITATION OF LIABILITY SHALL APPLY TO THE FULLEST EXTENT PERMITTED BY LAW.