



PRODUCT VALIDATION GRANT APPLICATION

APPLY NOW TO WIN a Product Validation Grant for your research.

Enhance your innovative research with Enzo's products.

Enzo scientists are on the lookout for innovative research projects that could benefit from our solutions.

[Check out Enzo's products and find out how to incorporate them into your workflow](#)

Winning entries will receive product kits free of charge and/or monetary grant to assist with the additional expenses to support the proposed project. To enter, fill out the application form and submit your entry.

FAQs

- **How do I apply?** Simply fill out the application form detailing your project and submit it along with your resume/CV to grants@enzo.com.
- **Who is eligible for this grant program?** The grant program is open to individuals working in academic research institutes. Other restrictions may apply.
- **What does the grant provide?** The principal investigator (PI) of the successful grant winner will receive product kits required to support applicant's research based on the project description (subject to submission of experimental plan). The grant may include funding for additional ancillary expenses.
- **When is the submission deadline?** All entries are reviewed on an ongoing basis.
- **What is the judging criteria?** Each entry will be judged by a team of scientists at Enzo on scientific merit, innovation, and fit with Enzo's technology.
- **What is required if I am awarded a grant?** The grant winner will be required to be available for participation in promotional activities in the first twelve months following announcement of the award. This could include conducting a webinar, assistance in write-up of a case study or application notes, or being a featured researcher in a spotlight/website/social media story. For more information, refer to the Terms & Conditions below. Additional travel award can be granted to present research at scientific conferences.



PRODUCT VALIDATION GRANT APPLICATION

Name: _____ **Degree/Title:** _____

Email Address: _____

Phone Number: _____

Institution: _____

Principal Investigator: _____

Job Title: _____

Department: _____

Address: _____

City: _____ **State:** _____ **Zip:** _____

Country: _____



*

Project Title: _____

Abstract:



Describe how you currently conduct DNA, RNA, and protein analysis. (Max. 200 characters with spaces)

*

Describe how you would incorporate Enzo's products into your project and workflow. Highlight how these assays would address current research challenges (max. 200 characters with spaces)

Please select your **research area(s)**:



- Bioprocess
- Cancer
- Cardiovascular
- Cell Signaling
- Cellular Biology
- Developmental Biology
- Drug Development
- Drug Discovery
- Epigenetics
- Genomics
- Immunology
- Immuno-Oncology
- Metabolism
- Molecular Biology
- Neuroscience
- Pathology
- Proteomics
- Proteostasis
- Spatial Biology
- Stem Cells
- Structural Biology
- Targeted Protein Degradation
- Toxicology
- Vaccine Development

SUBMIT APPLICATION TO: grants@enzo.com.

ENZO PRODUCT VALIDATION GRANT APPLICATION CHECKLIST

- Complete Product Validation Grant application
- 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 & Conditions

By submitting your entry, you agree to these terms for yourself, on behalf of 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.

Grant terms:

1. Only one entry per individual. Once entry is submitted, you may not amend it. Completed application and supporting documents must be submitted to qualify.
2. All qualified entries will be judged by a team of Enzo scientists for appropriateness and completeness in response. Each entry will be judged based on our view of how well the entry reflects scientific merit, innovation, and fit with Enzo's technology.
3. Successful grant applicants will receive product kits, controls, and ancillary products free of charge to support their project based on the project proposal description and subject to approval of the experimental plan by Enzo.
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.**

I understand and accept the terms and conditions.

By checking this box, you acknowledge that you read, understand, and fully accept the terms and conditions on this application. 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 on this application. To the extent you translated the aforementioned terms and conditions, you fully assume all risks associated with translating these terms and conditions.