

PBL Assay Science

Interferon Research Grant Application Form

Thank you for your interest in the PBL Assay Science Interferon Research Grant. One winner will receive \$10,000 in reagent credits for eligible PBL products. Credits may be used over multiple orders; shipping costs count toward the total. All research data generated remains the property of the recipient and institution. Completed applications should be submitted to: grant@pblassaysci.com

Applicant Information

Full Name:

Position/Title:

Institution Name:

Department / Lab:

Institution Address:

City:

State:

ZIP Code:

Email:

Phone:

Project Details

Project Title:

Project Summary (max 500 words):

Research Objectives:

Scientific Impact:

Innovation:

Workflow (Use of PBL reagents):



Budget & Resource Plan:



Agreements:

I agree to the Terms & Conditions.

I acknowledge all research data belong to myself and my institution.

I agree to participate in promotional activities if selected.

Upload up to two supporting documents (PDF preferred, max 10 MB each).

Submit completed application to: grant@pblsaysci.com