



**2023 IMMUNOLOGY2023™  
Annual Meeting Travel Award  
Application**

## Introduction

GeneCopoeia is pleased to announce the opening of a competition for an award for travel to and from the American Association of Immunologists (AAI) IMMUNOLOGY2023™ annual meeting, which will take place May 11-15, 2023 at the Walter E. Washington Convention Center in Washington, D.C. GeneCopoeia will be an exhibitor at the meeting to support our [OmicsArray™ line of antigen microarray products and services](#), and will provide the winner of the competition up to \$1,000.00 toward the travel expenses to and from the meeting.

Applicants must submit a description of their research, and how attending IMMUNOLOGY2023™ will benefit their research and contribute significantly to a field of immunology. Applications will be judged by a panel of GeneCopoeia scientists. The winning entry will be announced in early 2023.

To apply, please review the application eligibility requirements and the instructions, and complete the required information on the following pages.

## Eligibility requirements

To apply for the GeneCopoeia IMMUNOLOGY2023™ annual meeting travel award, applicants must meet the following eligibility requirements:

- At the time of the meeting, applications must be either graduate students enrolled at an accredited college or university, or b) employed as postdoctoral fellows.
- Applicants must be in an academic, government, or non-profit institution at the time of the meeting.
- Applicant's institution at the time of the meeting must be located within either the United States of America or Canada.

## Application period

- Application submissions begin Monday, August 1, 2022.
- The competition will close at 11:59 PM Eastern Time on Friday, October 28, 2022 or until 50 applications have been received, whichever comes first.

## Required documents

- This application.
- Curriculum vitae (CV) or résumé.
- Proof of enrollment as a Master's or Doctoral student at an accredited college or University within the United States of America or Canada between May 11-15 2023, OR employment as a postdoctoral fellow at an academic, government, or non-profit institution between May 11-15, 2023.

## Award delivery

The winning applicant will be reimbursed up to \$1,000.00 for expenses used for travel to and from the IMMUNOLOGY2023™ annual meeting May 11-15, 2023 in Washington, D.C. Submission of travel receipts to GeneCopoeia will be required.

**Instructions** (**Note:** Use of [Adobe Acrobat® Reader DC](#) is strongly recommended).

1. Fill out all fields in the “Personal information” section.
2. Provide a description of your research and how your participation in the IMMUNOLOGY2023™ annual meeting from May 11-15, 2023 in Washington, D.C. will benefit your research and make significant contributions to a field of immunology. Descriptions may not exceed 6,600 characters (approximately 1,000 words, including spaces, punctuation marks, and line breaks). Figures, tables, images, and/or data may be sent as separate attachments.
3. Save the form as a PDF.
4. Email the saved form, all required documents, and figures, tables, images, and data to [inquiry@geneCopoeia.com](mailto:inquiry@geneCopoeia.com). **Note:** Incomplete applications may lead to disqualification.

### Personal information (*all fields are required*)

Name of applicant: \_\_\_\_\_

Status of applicant (graduate student or postdoctoral fellow): \_\_\_\_\_

Applicant's institution: \_\_\_\_\_

Institution street address: \_\_\_\_\_

Institution city: \_\_\_\_\_

Institution state: \_\_\_\_

Institution ZIP code: \_\_\_\_\_

Institution email: \_\_\_\_\_

Institution phone number: \_\_\_\_\_

Name of Principal Investigator: \_\_\_\_\_

Email address of Principal Investigator: \_\_\_\_\_

**Description of research** (*required*)

## About GeneCopoeia

GeneCopoeia, Inc. a U.S.-based functional genomics company founded in 1999, develops cutting edge products and services for genomics, proteomics, and molecular and cellular biology. With emphasis on innovative R&D and superior product quality, GeneCopoeia serves the needs of biomedical scientists in academic research, and R&D of pharmaceutical and diagnostic products in their pursuit of discoveries and innovations.

Through the development, production, and marketing of the [OmicsArray™ line of antigen protein microarray products and services](#), GeneCopoeia is an industry leader in providing tools for autoimmune disease research. The progenitors of these arrays have been used for research by clinical and basic research laboratories that has led to dozens of high-impact, peer-reviewed publications.