



2024 GenScript Life Science

# Research Grant Program

Grant Application Form

# 2024 GenScript Life Science Research Grant Program Grant Application Form

## General information

Introducing the 2024 GenScript Life Science Research Grant Program.

Recognizing the pivotal role that research and innovation play in advancing the next generation of therapies and vaccines, GenScript is proud to announce the launch of the GenScript Life Science Research Grant Program—a new initiative dedicated to supporting breakthroughs in life science research areas including but not limited to gene and cell therapy development, antibody drug discovery, vaccine development, diagnostics, and plant sciences. This program is designed to empower researchers by providing grant funding specifically earmarked for purchasing GenScript reagents and services.

At GenScript, we understand the challenges and importance of advancing scientific knowledge of and developing treatments for life-altering diseases. In alignment with our commitment to fostering excellence in life sciences research, the GenScript Life Science Research Grant Program aims to catalyze innovation by offering financial support to researchers engaged in projects that can contribute significantly to critical areas of study.

Applications for this grant will be reviewed by a diverse committee comprised of GenScript's internal scientific leadership. Each proposal will be reviewed and ranked based on scientific merit, potential for impact, and strategic alignment with GenScript's primary areas of scientific interest.

The application process for the GenScript Life Science Research Grant Program will be transparent, efficient, and managed by a dedicated research grant team. Detailed information, including official rules, application procedures, documentation requirements, review processes, and decision notifications, will be available on our company website, [GenScript.com](https://www.genscript.com). GenScript is excited to contribute to advancing life sciences research by awarding grants at four pre-determined intervals throughout the year, fostering collaboration and excellence in research. The deadlines for each submission are as follows:

Deadline 1: 3/15/2024

Deadline 3: 9/15/2024

Deadline 2: 6/15/2024

Deadline 4: 11/15/2024

We expect to announce grant recipients within six (6) weeks of each submission deadline.

We look forward to receiving innovative proposals from researchers dedicated to making a positive impact in life science research areas including but not limited to cell and gene therapy, antibody drug discovery, vaccine development, diagnostics, or plant sciences. Together, let's

---

advance the boundaries of scientific knowledge and bring transformative solutions to the global life science, biotech, and medical communities.

## Application Instructions:

Grant applications for the 2024 GenScript Life Science Research Grant Program must be submitted electronically in English using the portal on the web page: <https://www.genscript.com/grantprogram.html>. Alternatively, completed application forms may be submitted to [LSRgrant@genscript.com](mailto:LSRgrant@genscript.com). Proposed projects should commence within six (6) months of receiving grant funding from GenScript, and research activities should generally take place over one year or less. Please note any planned deviation(s) from these guidelines in the application if applicable.

## Personal Information:

1. **Full Name:**
2. **Title/Position:**
3. **Institutional Affiliation:**
4. **Contact Information (Email and Phone):**
5. **Research Team Members:**
  - List names, titles, and roles of all team members involved in the project.

## Project Information:

6. **Project Title:**
7. **Research Area:**
  - Gene and Cell Therapy
  - Antibody Drug Discovery
  - Vaccine Development
  - Diagnostics
  - Plant Sciences
  - Others





---

#### Notice on Conflicts of Interest:

To ensure an unbiased and fair evaluation process, applicants are required to disclose any potential conflicts of interest that may compromise the impartiality of the review. Conflicts of interest may include, but are not limited to, financial relationships with GenScript, familial or professional relationships with members of the review committee, or any other affiliations that could influence the evaluation process. In the event of a conflict, GenScript reserves the right to take appropriate measures, including the disqualification of a proposal or the recusal of conflicted individuals from the evaluation process. GenScript is committed to upholding the highest standards of transparency, fairness, and confidentiality throughout the application and evaluation process for the GenScript Life Science Research Grant Program.

#### **14. Rights and Consent for Use of Participant Information and Project Data:**

To finalize award distribution, GenScript's Life Science Research Grant team will present a "Collaboration Agreement" for review to the grant recipients. The agreement aims to ensure that GenScript is publicly recognized in diverse forums for its contribution through the grant program. GenScript will not request that applicants share intellectual property (IP) rights to their research, findings, or products through this agreement. The awardee would own all rights in performing Services under the LSR Grant Program.

As a condition of participation, and in consideration for the opportunity to apply for the GenScript Life Science Research Grant Program, applicants agree to grant GenScript and its affiliates, and subsidiaries an irrevocable, perpetual, non-exclusive, and royalty-free right to use the participant's name, image, and likeness for any purpose and in all media, whether now known or hereinafter created, throughout the world without the expectation of compensation. Additionally, participants provide consent for GenScript and its affiliates to publish news releases regarding the GenScript Life Science Research Grant Program, grant recipients, and awarded research project titles on their websites and/or social media pages for promotional marketing purposes indefinitely. Participants further consent to being contacted by GenScript or its representatives for follow-up marketing activities related to the program. GenScript ultimately reserves the right to use any data generated from awarded research projects in marketing activities. However, GenScript will work with award recipients to determine the timing and content of such marketing activities.

---

## Supporting Documents (optional):

### 15. Attach any of the following documents:

- Research Proposal (PDF)
- Curriculum Vitae (CV) for each team member (PDF)
- Itemized Budget Request (PDF)
- Any additional relevant documents

## Declaration:

I hereby declare that the information provided in this application is true and accurate to the best of my knowledge. I understand that any misrepresentation may result in disqualification from grant consideration.

**Applicant Signature:** \_\_\_\_\_

**Date:** \_\_\_\_\_