



2025 GenScript Life Science  
**Research Grant Program**

Grant Application Form

# 2025 GenScript Life Science Research Grant Program Grant Application Form

## General information

Introducing the 2025 GenScript Life Science Research Grant Program.

Recognizing the vital role of research and innovation in driving the next generation of therapies and vaccines, GenScript proudly continues the GenScript Life Science Research Grant Program. This initiative is designed to empower researchers by offering grant funding specifically for purchasing GenScript reagents and services. Researchers from various life science fields are encouraged to apply and take advantage of this unique opportunity to advance their work.

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 two pre-determined intervals throughout the year, fostering collaboration and excellence in research. The deadlines for each submission are as follows:

Deadline 1: 6/15/2025

Deadline 2: 11/15/2025

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, and AI-driven drug discovery.

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 2025 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.
6. **Please provide a headshot (png, jpeg, or similar format) with your application.**

## Project Information:

7. **Project Title:**
8. **Research Area:**

- ☐ Gene and Cell Therapy
- ☐ Antibody Drug Discovery
- ☐ AI-Driven Drug Discovery
- ☐ Vaccine Development
- ☐ Diagnostics
- ☐ Others

---

**9. Project Scope (max 250 words):**

- Briefly describe the goals and objectives of your research project.

**10. Research Significance and Key Innovations (max 500 words):**

- Explain the importance and potential impact of your research in the chosen field and any novel approaches, technologies, or methodologies that set your project apart.

**11. Timeline (max 300 words):**

- Outline the milestones and key activities with corresponding timelines for your research.

**Budget and Funding:**

**12. Budget Breakdown (max 500 words):**

- This grant is designed to provide funding explicitly designated for purchasing GenScript reagents and services. Funds are not available to cover indirect costs, such as salaries, to buy products from other companies, or any other expenses.
- Provide your budget request amount and an estimated list of GenScript reagents and services awarded grant funds would be used for.
- The maximum budget for each project is set at USD 100,000.00. Only a few applications will be considered for maximum funding.
- Various grant funding levels will be considered, including 5,000-10,000, 10,000-50,000, and 50,000-100,000 USD. Applicants may request a specific funding level; however, GenScript reserves all rights to make the final decision on the funding amount awarded as part of the LSRG program.

**Eligibility and Compliance:**

**13. Eligibility**

*Research Focus:*

Proposals must align with the focus areas of the program in life science research areas including but not limited to gene and cell therapy development, antibody drug discovery, vaccine development, AI-driven drug discovery, or diagnostics.

*Geographic Location:*

Open to researchers globally. There are no restrictions based on geographic location, and international applicants are encouraged to apply.

---

#### *Entity Type:*

Eligible entities include academic/research institutions, universities, biotechnology companies, and other organizations engaged in life sciences research.

Any local, state, provincial, federal, or other government or government agency, and officials or employees thereof, are not eligible to participate.

Healthcare organizations, including hospitals, healthcare professionals, and employees thereof, are not eligible.

#### *Legal Status:*

Must be legally recognized and authorized to conduct research in their respective jurisdictions.

#### *Affiliation Restrictions:*

Employees of GenScript or its affiliates, as well as their immediate family members and household members, are ineligible to apply.

#### *Collaborations:*

Each eligible entity may submit only one (1) application representing the sole efforts of that entity. Collaborations among two (2) or more entities will not be considered unless GenScript provides explicit written consent.

#### *Compliance:*

Applicants must comply with all applicable federal, state, and local laws and regulations.

#### *Void Where Prohibited:*

The program is open to participants worldwide, except where prohibited by law. Participation constitutes the applicant's full and unconditional agreement to the program's rules and decisions, which are final and binding in all matters related to the grant.

GenScript reserves the right to verify the eligibility of applicants and to disqualify any application that does not meet the specified criteria. The eligibility criteria aim to ensure a fair and open application process for researchers across diverse backgrounds and geographic locations.

#### *Terms and Conditions:*

GenScript reserves the right to make the final determination of funding for the Life Science Research Grant (LSRG) Program. The selection of individuals to receive LSRG awards, the total number of awards, and the amount of funding provided are solely determined by GenScript. Applicants may request a specific budget based on the proposal submitted during the application process; however, GenScript retains all rights to decide on the funding level and the extent of support provided to each awardee toward services and reagents.

---

#### **14. Confidentiality and Conflicts of Interest:**

##### **Confidentiality Terms:**

GenScript recognizes the sensitive nature of research proposals and respects the proprietary information submitted by applicants. Therefore, all details provided in the grant applications, including methodologies, data, and preliminary findings, will be treated as confidential.

##### **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.

#### **15. 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):

### 16. 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:** \_\_\_\_\_