

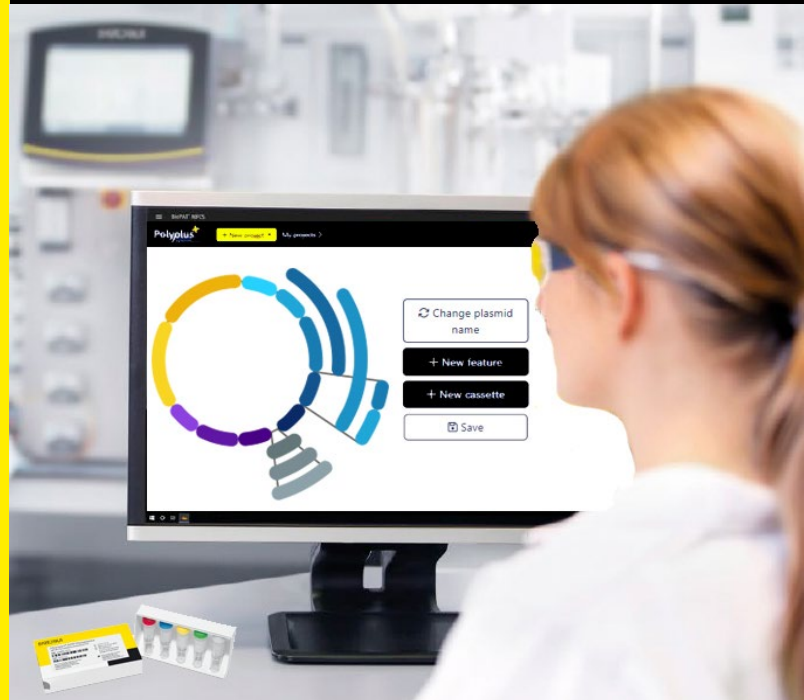
Easy Plasmid Service

Plasmid Designed to Cater To Your Specific Requirements, Ensuring a Seamless And Cost-Effective Experience

Service Information

Our plasmid platform offers a range of easy-to-use plasmids that can provide you with several benefits for your scientific projects. Based on e-Zyvec® technology, we create a library of plasmid template to fit your needs. Use our online platform to choose the plasmid that best suits you, modify your expression cassette, and submit your request.

With the help of our experts, get a review of your plasmid before placing an order, and then receive your ready-to-use plasmids in the desired quantity for your project.



Features and Benefits

- **Cost-Effective:** Reduce your project costs by using easy plasmids (from 249 €)
- **Scalable:** Access highly complex plasmids with our tailor-made service, up to GMP grade
- **Quality Assurance:** Plasmid scientist experts and whole plasmid sequencing included
- **Time-Saving:** Ready-to-use discovery grade plasmids within 5 weeks from ordering

Introduction

Relevant Applications

- Fundamental research
- Cell and gene therapy: AAV and LV transfection

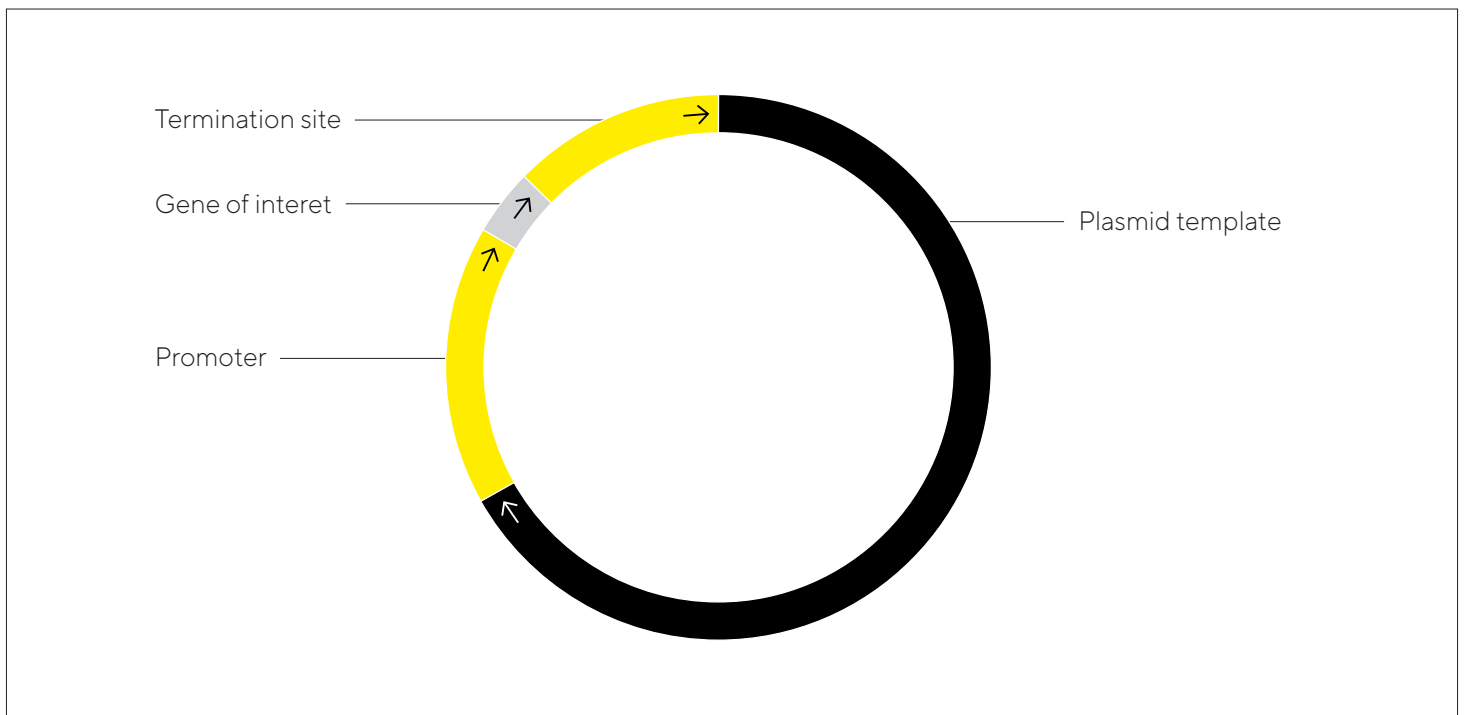
Quality Grade Provided

Easy plasmid service provides plasmid amplification up to several mg of plasmids as discovery grade (Miniprep and whole plasmids sequencing as QC are included).

Higher quantity and | or quality (up to GMP grade) can be performed through our manufacturing services.

Easy Plasmid Service Workflow

- 1. Access:** Log in to our online design platform and start a new plasmid project.
- 2. Select:** Choose an easy plasmid template that meets your requirements (if any plasmid architecture fits, create a tailor-made one through our plasmids engineering service, nothing is restricted).
- 3. Design:** Modify the expression cassette by selecting a promoter and inserting your gene of interest.
- 4. Review:** Submit your design for a complimentary expert plasmid review.
- 5. Confirm:** Approve the final design and specify the required plasmid quantity.
- 6. Production:** Keep updated on the production progress of your plasmids through our online platform.
- 7. Quality Control:** Validation of your plasmid thanks to different QCs as whole plasmids sequencing.
- 8. Move Science Forward:** Finalize your project with our ready-to-use plasmids.



Ordering Information

Easy plasmids have been designed to make the process of plasmids design and production, easier and cheaper (from 249 €).

To proceed with the order, create an account at www.polyplus-sartorius.com/register, then access the design platform. Create your project and choose your easy plasmids, then submit your plasmids project. It will be reviewed by our experts, who will return to you with their suggestions. Then a quote will be issued.

Product Name	Product Description	Order Number
Plasmid pTransgene AAV	Plasmid Transgene to be co-transfected into packaging cells along with the pHelper plasmid (e.g. pPLUS [®] AAV-Helper), the pRepCap plasmid, and transfection reagents (e.g. FectoVIR [®] -AAV), enabling the efficient production of viral particles to further transduce targeted cells tissues to express the gene of interest.	101000221
Plasmid pTransgene LV	Plasmid Transgene plasmid to be co-transfected into packaging cells along with packaging lentiviral plasmids encoding the Gag, Pol, Rev, VSV-G viral components, and with the transfection reagents (e.g. FectoVIR [®] -LV).	101000220
Plasmid shRNA Expression	The easy shRNA plasmid to base-pairs with a complementary target messenger RNA (mRNA), leading to its degradation.	101000219
Plasmid Protein Expression	The easy protein expression plasmid engineered by Polyplus carry the genetic features required by mammalian systems to synthesize and express the desired protein.	101000218

End Product Information

We deliver discovery grade plasmids as circular, double stranded DNA molecules, in a screw-cap plastic tube, re-suspended in 1 × AE buffer (miniprep.) or 1 × TE buffer (midi-, maxi-, giga- preparation.), at room temperature. Please note that we do not provide bacterial stock.

For research-, high quality- and GMP-grade, please refer to our plasmid manufacturing service.

Quality Controls

QCs are limited to basic quality and identity tests:

- DNA quantification
- A260/280 and A260/230 measurements
- Restriction digest profile (gel or capillary electrophoresis)
- Whole plasmid sequencing including mycoplasma and gDNA detection (NGS)
- Sanger sequencing (if necessary)

QC results will be documented in the Product Delivery Form. Endotoxins levels are not measured for discovery grade plasmids.

Storage

Upon receipt, briefly spin down the tubes containing your engineered plasmids (bench centrifuge) and store them at:

- 4 °C for short storage (<1 month)
- -20 °C for long storage

We strongly recommend users aliquot discovery grade plasmids for long storage and to avoid freeze-thaw cycles. Duration of storage is given here as an indication. Always verify the integrity of plasmids stored for extended periods by gel electrophoresis (denaturation) or sequencing (sequence integrity).

Intellectual Property

Unless otherwise stated in the specifications of you project, every engineered plasmid provided on a service based contract, for Research Application, are IP free from Sartorius, as stipulated in our T&C. Other applications may require licensing agreement with Sartorius or tiers entity.

Germany

Sartorius Stedim Biotech GmbH
August-Spindler-Strasse 11
37079 Goettingen
Phone +49 551 308 0

USA

Sartorius Stedim North America Inc.
565 Johnson Avenue
Bohemia, NY 11716
Toll-Free +1 800 368 7178

France

Polyplus®—Now part of Sartorius
75 Rue Marguerite Perey
67400 Illkirch
Phone +33 390 406 180



For more information, visit

sartorius.com/transfection-reagents-plasmids