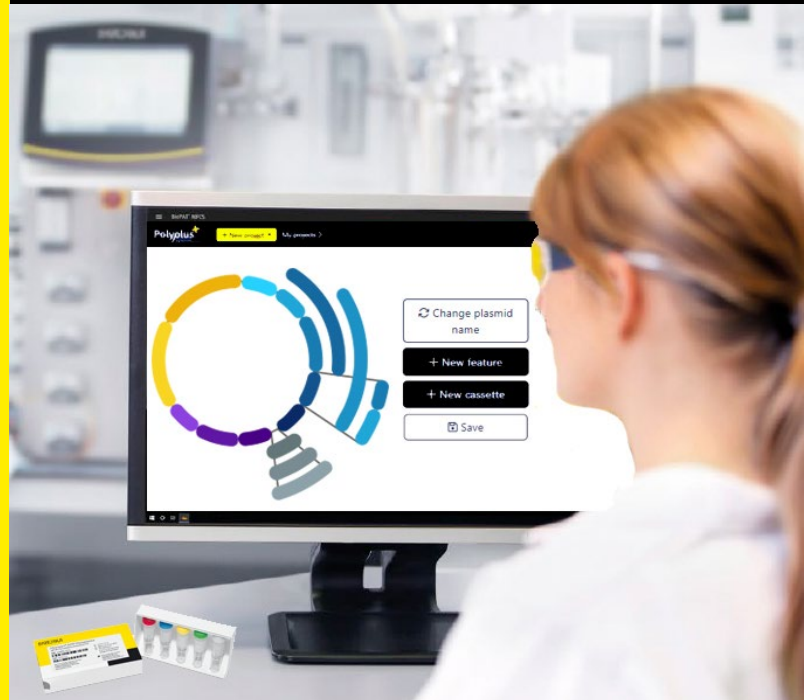


Tailor-Made Plasmid Service

Engineering and Production Services for Plasmids From Discovery to GMP-Grade



Service Information

Our unique and innovative DNA assembly technology, e-Zyvec®, utilizes linear DNA bricks to design and assemble tailor-made plasmids. This allows for fully flexible and accurate plasmid design, enabling the generation of any plasmid, from simple to highly sophisticated, without restrictions.

Using our online plasmid platform, you can design your plasmid, share it with your team, and consult with our experts who can help identify your needs from the start of your project. You will receive ready-to-use, tailor-made plasmids that meet your project's expectations. Additionally, the full length of your plasmid will be validated by next-generation sequencing as a quality control measure.

Features and Benefits

- **Innovative Technology (e-Zyvec®):** Seamless DNA assembly to produce even highly complex plasmids
- **De-Risking:** Access to dedicated plasmid experts and an online design platform
- **Reliable Plasmid Sequence:** Full-length plasmid sequencing included
- **Time-Saving:** Ready-to-use plasmids within three weeks of ordering

Introduction

Relevant Applications

- Fundamental research
- Cell and gene therapy
- mRNA and DNA therapeutics
- Protein and antibody engineering

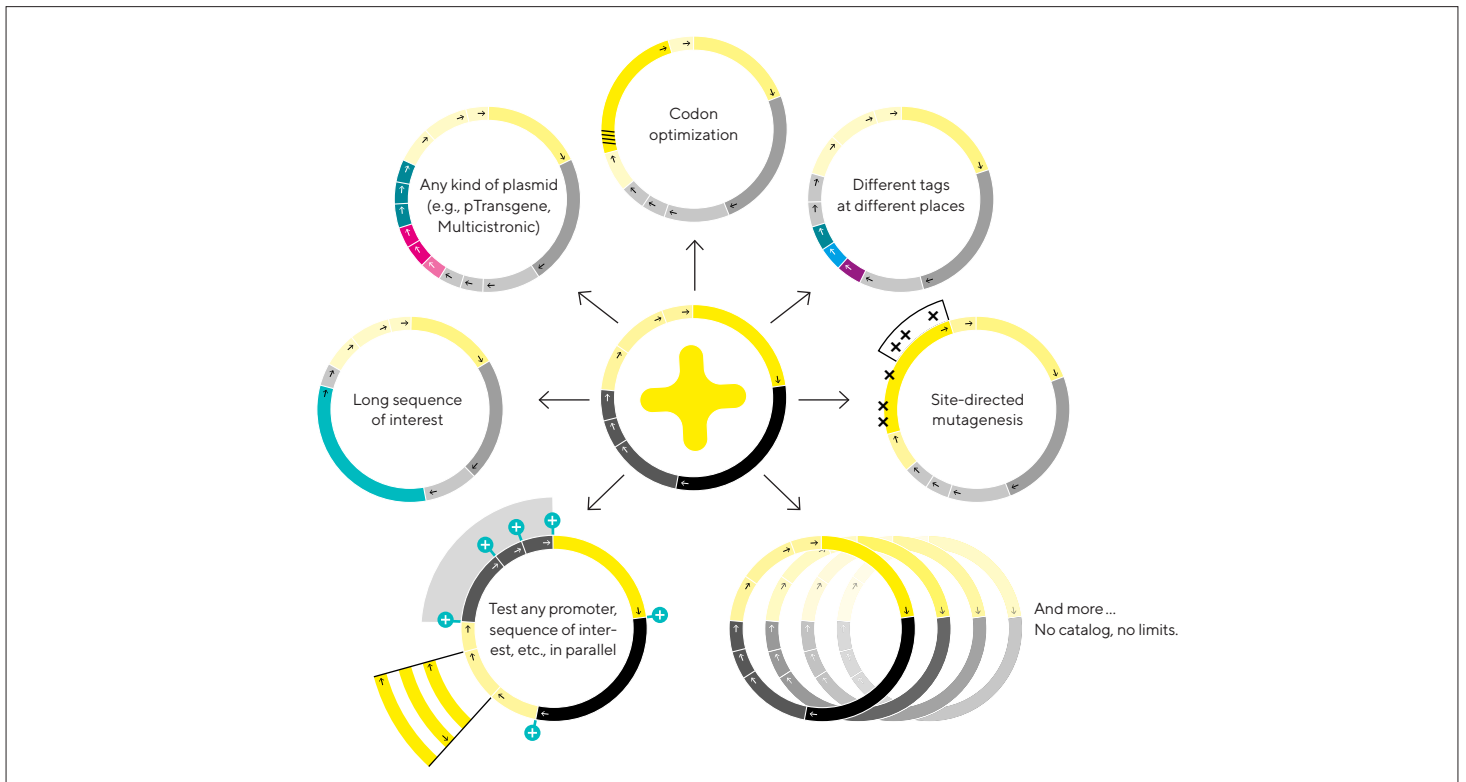
Quality Grade Provided

Our tailor-made plasmid service provides plasmid amplification up to several milligrams at discovery grade (miniprep and whole-plasmid sequencing are included as quality control).

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

Tailor-Made Plasmid Service Workflow

- 1. Access:** Log in to our online design platform.
- 2. Initiate:** Start a new plasmid project.
- 3. Select:** Choose to design a plasmid from scratch or use a suggested template.
- 4. Design:** Start to assemble a plasmid from scratch or use the suggested template. Add and move any sequence from our database. Everything is flexible, nothing is restricted.
- 5. Review:** Submit your design for a complimentary expert plasmid evaluation.
- 6. Confirm:** Approve the final design and specify the required plasmid quantity.
- 7. Track:** Stay updated on the production progress of your plasmids through our online platform.
- 8. Quality Control:** Your plasmid is validated thanks to different quality control measures, including whole-plasmid sequencing.
- 9. Move Science Forward:** Finalize your project with our ready-to-use plasmids.



Ordering Information

To proceed with the order, register at www.polyplus-sartorius.com/register, create a project, and start designing your plasmid with our easy-to-use platform.

After designing your plasmid on our online platform, you can get in touch with our plasmid scientist to review your plasmids and get advice! Once you validate the design, our experts and our proprietary software will engineer your plasmid to maximize your chances of success (99% success rate).

For tailor-made plasmid projects, desired sequences can typically be generated using DNA synthesis (outsourced). However, some projects may require you to send us DNA material (plasmid templates or purified genomic DNA) that we can use to overcome technical complexity, ensure technical continuity, or reduce cost. In this case, please follow these instructions:

- Use a 1.5 mL plastic tube, hermetically closed (screw cap or parafilm)
- Label the tube clearly
- Minimal volume: 15 μ L
- Minimal concentration: 100 ng/ μ L
- Room temperature, please avoid dry ice

End Product Information

We deliver discovery grade plasmids as circular, double-stranded DNA molecules, in a screw-cap plastic tube, re-suspended in 1 \times AE buffer (miniprep) or 1 \times 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

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 genomic DNA detection (by next-generation sequencing)
- Sanger sequencing (if necessary)

Quality control results will be documented in the Product Delivery Form. Endotoxin 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 $^{\circ}$ C for short storage (<1 month)
- -20 $^{\circ}$ 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 your project, every engineered plasmid provided on a service-based contract for research applications is IP-free from Sartorius, as stipulated in our T&C. Other applications may require a 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