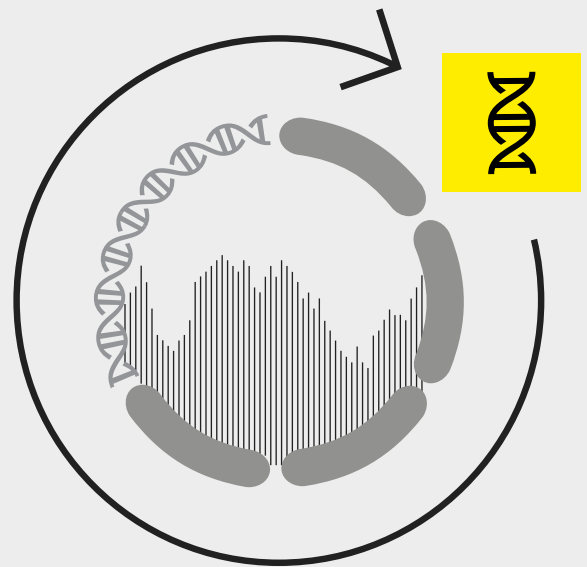


Whole Plasmid Sequencing Service

Full-Length Sequencing of Multiple DNA Plasmids



Service Information

Our state-of-the-art Whole Plasmid Sequencing Service provides high-quality and accurate sequencing of any plasmid, irrespective of size and complexity.

The service delivers rapid results (from three working days upon receipt of plasmids). Beyond the data, our personalized consultations and competitive pricing make plasmid sequencing accessible to all researchers, with the possibility to sequence a single plasmid sample, or up to 150 samples in one run.

Features and Benefits

- **Time-Saving:** Receive your whole plasmid sequencing results in as little as three working days.
- **Personalized Reports:** Full-length sequencing of predictive (sequence matching) or unknown plasmids (de novo analysis)
- **Expertise:** In-house bioinformatical pipeline allowing quantification and identification of contaminant sequences

Introduction

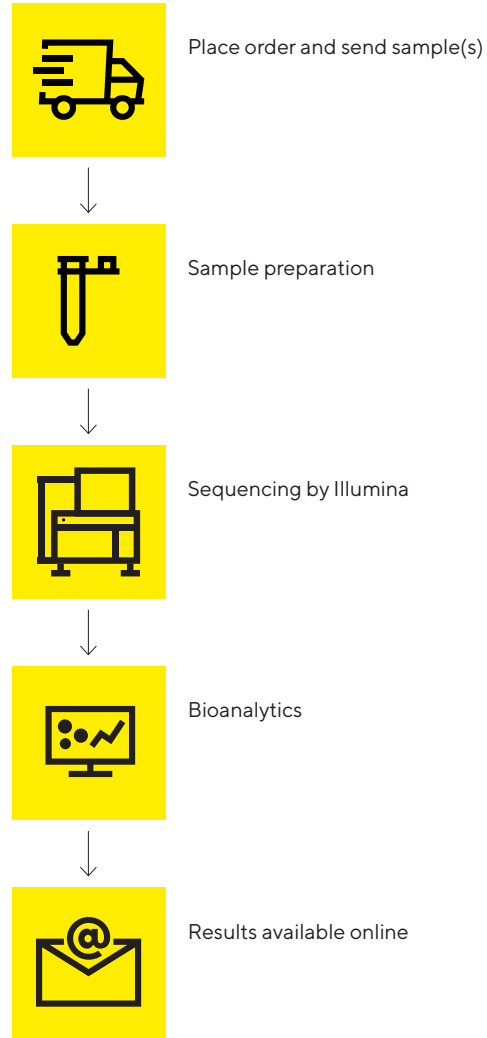
Our Expertise

We have spent years refining our understanding of plasmid structures, sequences, and variations. This expertise enables us to deliver precise insights, including de novo analysis, that advance your research and applications. Our capabilities extend beyond data provision; we also partner with you to navigate the complexities of plasmid-related challenges. Thanks to our service, you get more than a sequencing service—you access a depth of plasmid expertise designed to elevate your scientific journey.

Delivery Time

Getting your sequencing results is now faster than ever. Ordering sequencing for 12 or more plasmids guarantees the availability of results within three working days of sample receipt, ensuring rapid progress for your projects. Alternatively, our standard service delivers comprehensive results as early as two weeks following sample receipt. This fast turnaround time allows you to quickly obtain the sequence information you need to move forward with your research projects.

Plasmid Sequencing Workflow



Ordering Information

Our plasmid sequencing services are dedicated to whole plasmid sequencing. Therefore, the starting material should be purified double-stranded DNA plasmids. We offer next-generation sequencing for even a single sample, with no upper limit on the number of samples – we can sequence up to 150 samples simultaneously. Please note that we do not perform whole genome sequencing of any organism.

After sending an order through our online platform, you can send your samples following the guidelines below:

- 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

Samples should be sent to:

Polyplus-Transfection
Production Team
80 rue du Dr Yersin
59120 Loos
FRANCE

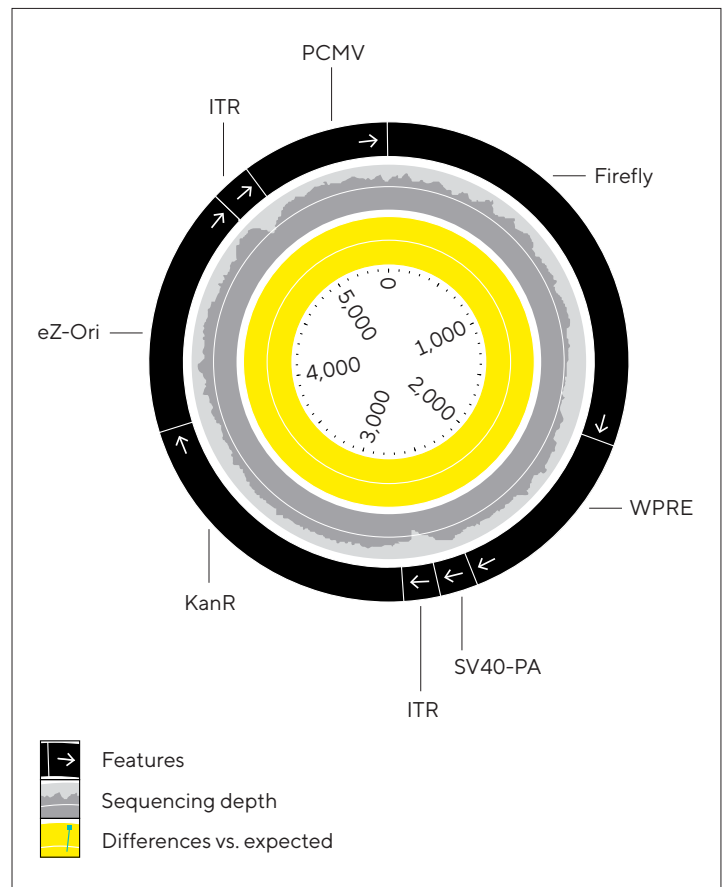
When known, please enter the sequences of reference of your samples on our platform.

Deliverables

After sequencing your sample, we will provide the full-length sequence of your plasmid (.fasta file) as well as a sequencing report (.pdf file) containing:

- A visual representation of the sequencing results with the coverage and the differences observed with the sequence of reference, if there is any
- The positions and the nature of the variations (insertion, deletion, substitution) observed in comparison with the sequence of reference
- The percentage of contaminations contained in your sample, such as bacterial gDNA

The sequencing report and the full-length sequence of your plasmids are available on your account.



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