

Laying the Foundation for Scientific Progress and Innovation *together*





Product Portfolio

DNA & RNA Analysis

- ▶ Cloning, Enzymes, and DNA Markers
pGEM-T®
- ▶ Transfection and Epigenetics
FuGENE® | ViaFect™
- ▶ DNA and RNA Purification
Maxwell® | PureYield™ | ReliaPrep™
- ▶ DNA and RNA Quantification
QuantiFluor® Dyes
- ▶ DNA Amplification, PCR and qPCR
GoTaq®
- ▶ Reverse Transcription
GoScript™
- ▶ RNase Inhibitors
RNasin®
- ▶ NGS Sample Preparation
ProNex®
- ▶ Sanger Sequencing
Spectrum Compact | ProDye®
- ▶ STR Analysis
PowerPlex® | GenePrint®

Cell Biology

- ▶ Gene and Protein Reporters
NanoLuc® | Ultra-Glo™ | Nano-Glo®
- ▶ Viability, Cytotoxicity and Apoptosis
CellTiter-Glo® | CellTox™
- ▶ Cell Metabolism
Glucose-Glo™
- ▶ Oxidative Stress
ROS-Glo™
- ▶ 3D Culture Analysis
Caspase-Glo® 3/7 3D
- ▶ Real-time Analysis
RealTime-Glo™
- ▶ Signal Transduction
RapidResponse™
- ▶ Cellular Imaging
HaloTag®
- ▶ Immunoassays
Lumit™

Protein Analysis

- ▶ Protein Expression
TNT® | KRX Cells
- ▶ Protein Quantification
Lumit™ | HiBiT
- ▶ Protein Degradation
NanoBRET™ | HiBiT | HaloTag®
- ▶ Pull-down Assays
MagneGST™ | HaloTag®
- ▶ Protein Interaction
NanoBiT® | NanoBRET™ | Lumit™
- ▶ Protein Purification
MagneBeads® | HaloTag®
- ▶ Mass Spectrometry
Trypsin/Lys-C | Trypsin Gold | AccuMAP™

Instruments

- ▶ Luminometer, Fluorometer and Multimode-Reader
GloMax® Detection Systems
- ▶ Automated DNA and RNA Extraction
Maxwell® | Maxprep™
- ▶ DNA and RNA Concentration Measurement
Quantus™ Fluorometer | GloMax® Detection Systems
- ▶ Capillary Electrophoresis for Sanger Sequencing and Fragment Analysis
Spectrum CE Systems

Reporter Potency Bioassays

- ▶ Fc Effector Function
ADCC, ADCP
- ▶ Cytokines & Growth Factors
VEGF, IL-2, IL-6, IL-12 and many more
- ▶ Immune Checkpoints
PD-1/PD-L1, CTLA-4, LAG-3, TIGIT
- ▶ T Cell Activation
Bioassays & knock-out variants (TCRαβ-KO, CD4+, CD8+)
- ▶ Drug & Target:
Interaction and Activity
PROTACs | IMiDs | Molecular Glues & GPCRs | Kinases | PARPs | Ras



www.promega.com/products



Explore Promega

Driven by curiosity and inspired by research, we have been supporting your success in science for over 40 years.

Promega promotes biological research and development worldwide with innovative products, expert support, and versatile services. The product diversity includes proven reagents and instruments for gene, protein, and cell analysis. With a global workforce of nearly 2,000 employees, we proudly serve customers in over 100 countries. Working in partnership and supporting scientific questions is our highest priority. Together, we want to leverage the full potential of life sciences and contribute to an overall improvement of the quality of life.

Everything from a single source: Reagents – Instruments – Service

We want to make our customers' work as easy as possible by offering complementary products and services. Our offerings includes expert product support, in-house manufacturing of customized products, batch reservations, instrument service, and training programs.



Your Application – Our Expertise



Fundamental Research

- ▶ Cell and Molecular Biology
- ▶ Immunology
- ▶ Genetics and Genomics
- ▶ Proteomics
- ▶ Epigenetics
- ▶ Biochemistry
- ▶ Systems Biology
- ▶ Physiology



Genetic Identity

- ▶ Forensic Trace DNA Analysis
- ▶ Human Identification
- ▶ Relationship Testing
- ▶ Capillary Electrophoresis
- ▶ Cell Line Authentication



Animal Health

- ▶ Pathogen Detection
- ▶ Disease Diagnostics
- ▶ Vaccine Development and Production
- ▶ Genotyping

Pharmaceutical Research & Development

- ▶ Small Molecule Drug Discovery
- ▶ Therapeutic Antibodies
- ▶ Cell, Gene and Immunotherapy
- ▶ Vaccine Development
- ▶ Customized Screening, Potency, and MoA Analyses



Environmental & Food Analysis

- ▶ Pathogens in Water, Plant, and Environmental Samples
- ▶ Epidemiological Wastewater Monitoring
- ▶ Plant Genotyping
- ▶ Food Authentication
- ▶ Feed and Seed Analysis



Clinical Research & Diagnostics

- ▶ Molecular Diagnostics, also as IVDR Compliant Solution
- ▶ Molecular Oncology
- ▶ Microsatellite Instability Testing (MSI)
- ▶ Drug Development



www.promega.com/applications

Service For You

Your success drives our commitment to service.

Technical Service

Scientific support with product selection and experimental setup

Comprehensive troubleshooting

Reliable and fast support

Application-specific protocols

Training Events

Seminars, webinars, and workshops on new technologies, products and applications

Promega Academy: individual seminars for your institute – live or online

Advanced trainings on laboratory methods and other topics



Instrument Service

In-depth consultation and financing options

Instrument installation and on-site training

Service contracts for maintenance and warranty

Maintenance and repair at the instrument service center in Germany or at your laboratory

Automation Support

Solutions for specific automation problems

Implementation of Promega products on liquid handlers of well-known manufacturers

On-site service & hands-on support

Product Innovations

Access to the latest product developments prior to market entry

License-free commercial use of many products

Customized Products

Individual filling and packaging options

Customized product formulations

Large-scale production / Custom OEM

Assay & cell line development



Helix® – Your On-Site Stocking System

Storage of Promega products in the laboratory

Freezers and room-temperature cabinets with intelligent RFID-based billing technology

Resource saving through consolidated deliveries and efficient refilling plans

Delivery & Supply Chain Security

Short delivery time from the European central warehouse in Germany

99% of all products are available immediately

Consistent quality through certified logistics processes

Optimized raw material and inventory management as well as in-house production



Storage Space & Batch Reservation

Storage options for customized products

Reservation of specific batches



www.promega.com/support



Exclusive Instrument Service

We provide comprehensive instrument service:
from detailed consultation and demonstration to installation
and after-sales support.

Consulting & Instrument Demonstrations

Get detailed guidance on instrument's functions. We are happy to demonstrate all instrument functions to you either on site in your laboratory, online or in our training room in Walldorf, Germany.

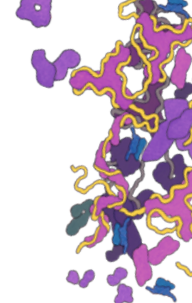
Purchase, Rent, or Financing

To facilitate access to an instrument of your choice, various purchase, rental, and financing models are available. Please contact us for consulting and support.

Delivery

Instrument delivery aligned to your working schedule.





You can rely on us:

We provide cost transparency and ongoing support throughout the entire instrument lifecycle.

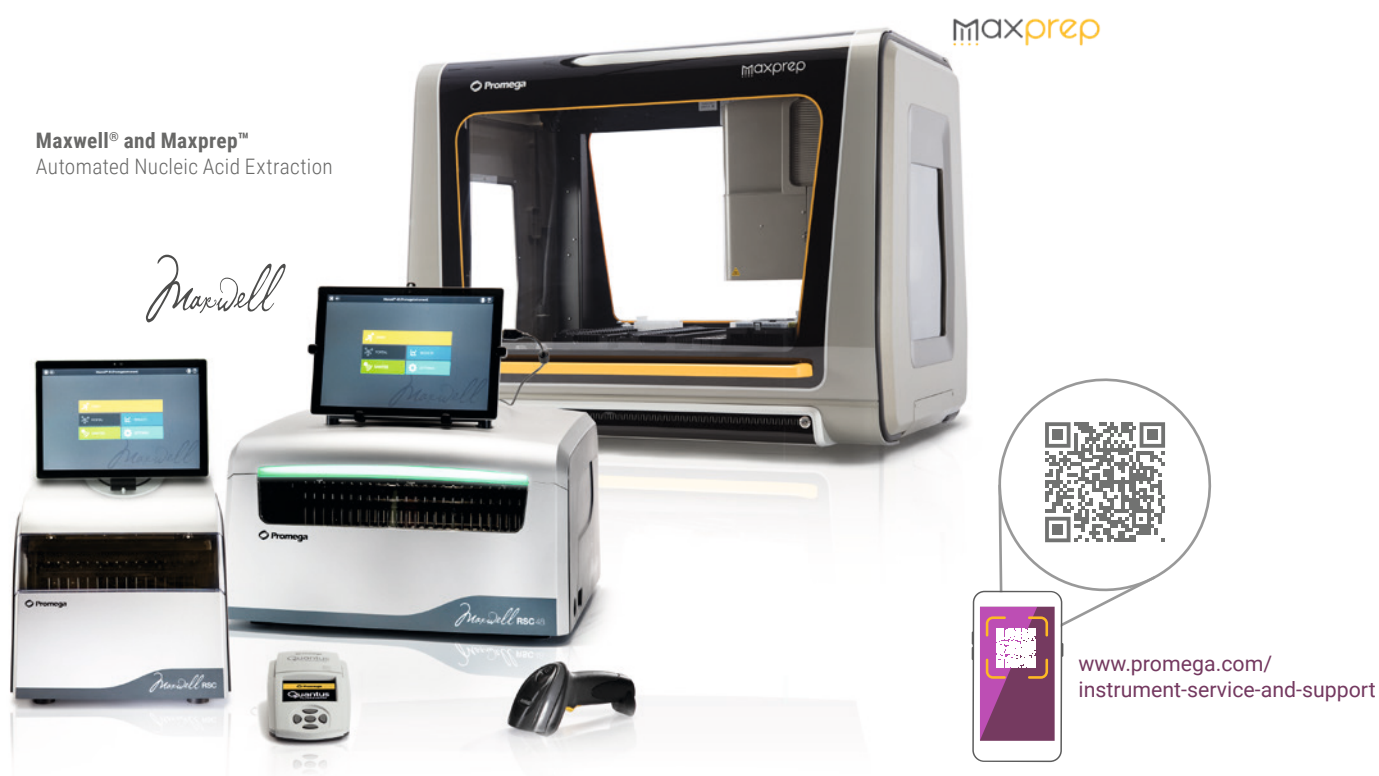
Installation & Training

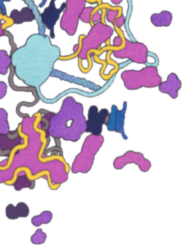
The instrument installation will be carried out in close coordination with your Promega Account Manager. You will be trained in the operation of the instrument by an application specialist according to your needs.

After-Sales Support

After successful installation of your instrument, our technical service will address your questions related to instrument operation and maintenance, reagents, new protocol design and implementation.

Instrument maintenance is provided by the service center in Germany. You will be notified in time about upcoming maintenances needed and expiring warranties. A variety of service contracts with different service options are available.





Luminescent Technologies

Our expertise for your applications – benefit from first-hand support with the development of bioluminescent test systems.

Our custom assay service offers you tailored assay development for your drug discovery, development, and QC workflows. This ranges from de novo assay design, modification of existing assays, assay optimization/qualification, to cell line manufacturing (e.g., using CRISPR) and drug profiling screening services.



- ▶ **Luminescent Reporters for Assay Development** based on HaloTag® labeling and NanoLuc® and Firefly luciferase technologies
- ▶ **Lumit™ Immunoassays** ready-to-use formats and options to build customized assays
- ▶ **NanoBiT® / NanoBRET™ Technologies** for protein:protein interaction assays in living cells. Suitable for drug screening of protein interaction modulators
- ▶ **NanoBRET™ Target Engagement Assays** to assess drug permeability and to characterize drug:target interactions
- ▶ **HiBiT Protein Tagging System** for the development of cell-based protein reporter assays that are also suitable for kinetic analyses in living cells. Protein detection is performed without antibodies

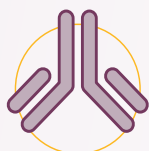


www.promega.com/tailored-solutions

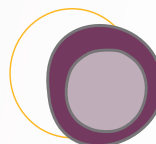


We are pleased to support you in implementing luminescent technologies in your research projects. Are you working on therapeutic antibodies, vaccines, cell or gene therapeutics? We will help you to develop the appropriate assays that reflect your drug's mode of action.

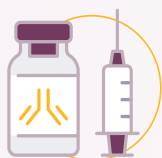
Therapeutic
Antibodies



T-cell Therapy



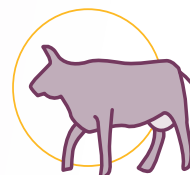
Vaccines & Antiviral
Antibodies



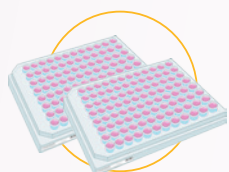
Our service for you:

- ▶ Potency assay development
- ▶ T-cell receptor discovery
- ▶ Bioassay qualification
- ▶ Drug profiling
- ▶ Cell production on request

Animal Health



Screening Service



Gene Therapy



Sustainability & Social Responsibility

Environmental Protection

Promega invests in numerous initiatives and projects to protect the environment. Examples include:

- ▶ Increased use of renewable energy to reduce consumption of energy from fossil fuels
- ▶ Optimized and recyclable product packaging for waste reduction
- ▶ Increased use of rainwater for water conservation
- ▶ Reduction of CO₂ emissions through the Helix on-site stocking system at customer's site, and much more

We are constantly working to improve our eco-balance. Our goal for 2030 is to reduce waste and water consumption by 30% and CO₂ emissions by 50% (based on fiscal 2019).



[www.promega.com/
corporate-responsibility-csr](http://www.promega.com/corporate-responsibility-csr)



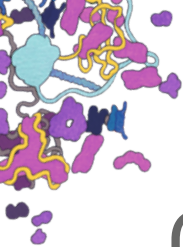
♥ Committed to Our Fellow Human Beings

We actively integrate the principles of sustainability throughout our company.


We are building and maintaining long-term relationships around the world by engaging local communities. Supporting education and science is a fundamental part of this.


These are some examples:

- ▶ Sponsoring young scientists to participate in the iGEM competition
- ▶ Coordinate events for employees to raise money or in-kind donations for local charities
- ▶ Support for an agricultural research project in Kenya
- ▶ Encourage minorities through equal opportunities in the hiring and promotion process



Contact

 Headquarters

 R&D and Production

 Branches

 Distributors

Europe

Germany/Austria

Tel +49 6227 6906 0
Fax +49 6227 6906 222
de_custserv@promega.com

Switzerland

Tel +44 878 90 00
Fax +44 878 90 10
ch_custserv@promega.com

France

Tel +33 0437 2250 00
Fax +33 0437 2250 10
contactfr@promega.com

Italy

Tel +39 0254 0501 94
Fax +39 0256 5616 45
customerservice.italia@promega.com

Poland

Tel +48 22 531 0667
Fax +48 22 531 0669
pl_custserv@promega.com

Central/Eastern Europe

Tel +49 6227 6906 0
Fax +49 6227 6906 222
de_ceesales@promega.com

Belgium/Luxembourg/Netherlands

Tel +31 71 532 42 44
Fax +31 71 532 49 07
benelux@promega.com

Spain

Tel +34 902 538 200
Fax +34 902 538 300
esp_custserv@promega.com

Denmark, Estonia, Finland, Iceland, Norway, Sweden

Tel +46 8 452 2450
Fax +46 8 452 2455
sweorder@promega.com

United Kingdom

Tel +44 23 8076 0225
Fax +44 23 8076 7014
ukcustserve@promega.com

North America

USA/Canada

Tel +1 608 274 4330
Fax +1 608 277 2516
custserv@promega.com

South America

Brazil

Tel +55 11 5090 3780
Fax +55 11 5096 3780
promega.brasil@promega.com





Asia

China

Tel +86 10 5825 6268
Fax +86 10 5825 6160
info@promega.com.cn

Japan

Tel +81 3 3669 7981
Fax +81 3 3669 7982
jpmktg@jp.promega.com

Korea

Tel +82 2 1588 3718
Fax +82 2 2638 5418
CustServiceKR@promega.com

Asia-Pacific region, Singapore

Tel +65 6513 3450
Fax +65 6773 5210
sg_custserv@promega.com

India

Tel +91 11 43005814/15/16/17
Fax +91 11 41035028
ind_custserv@promega.com

Australia

Australia

Tel +61 2 8338 3800
Fax +61 2 8338 3855
auscustserv@promega.com

More information about Promega
branches and distributors:

**[www.promega.com/support/
contact-promega](http://www.promega.com/support/contact-promega)**



