



Lab Automation 101 from Concept to Reality

Expert Implementation & HT Solutions



Overview

- » **Is your lab's efficiency suffering due to high workloads? Need higher throughput to keep up?**
- » **Facing challenges meeting regulatory standards while ensuring consistent, reproducible data?**

Automating your lab unlocks valuable benefits—saving time, increasing throughput, and improving quality. But getting there presents various challenges, from selecting the right platform and chemistry to integrating everything seamlessly. The abundance of available options can be overwhelming.

However, you don't have to face the complexities of automation implementation alone. When you partner with Promega, you gain access to our team of Field Support Scientists and Field Applications Scientists. With a combined experience of over 125 years in automation chemistries and platforms, our expert team is here to guide you efficiently from concept to a fully functional workflow.

And our support doesn't stop there—we remain accessible whenever you need assistance, offering expert guidance in both biology and automation. When you choose Promega, you're investing in a partnership dedicated to guaranteeing your success in achieving a seamless and successful automation transition. Trust us to be your dependable and capable partner every step of the way.

- » Excited to start with automation but uncertain about where to initiate your lab automation journey?
- » Overwhelmed by the multitude of options available?
- » Unsure of the best path forward?



Reach out for expert support and tailored recommendations during every step of your automation journey.

Learn how our FSS team helped to streamline high-throughput nucleic acid extraction, enabling efficient SARS-CoV-2 testing alongside existing workflows.



Inquire about Expert Automation Support

www.promega.com/labautomation



Benefits

Increased Efficiency and Throughput:

Accelerate experimentation and analysis, handling larger sample volumes in less time.

Enhanced Accuracy and Precision:

Minimize human error for consistently accurate and precise results.

Time and Resource Savings:

Free up your researchers' time and reduce labor costs with automated, repetitive task handling.

Cost-Effective Operations:

Initial investment yields long-term savings through minimized labor costs and error prevention.

Improved Data Quality and Documentation:

Generate standardized, well-documented data, ensuring compliance and easy reproducibility.

Expanded Experimental Capabilities:

Execute intricate experiments impractical or impossible manually, unlocking new research avenues.

Optimized Workflow Integration and Adaptability:

Create a seamless, interconnected workflow with integrated lab automation processes. Easily adapt or expand automated systems to meet evolving research needs and workloads.



Applications

Nucleic Acid Analysis



- ▶ gDNA purification
- ▶ RNA purification
- ▶ Plasmid extraction
- ▶ ccfDNA purification from plasma
- ▶ STR workflows
- ▶ DNA & RNA Clean-up
- ▶ Preparation of NGS Libraries

Protein Analysis



- ▶ Purification of antibodies
- ▶ Protein sample preparation for mass spectrometry
- ▶ Purification of polyhistidine proteins
- ▶ Purification of biotinylated antibodies and proteins

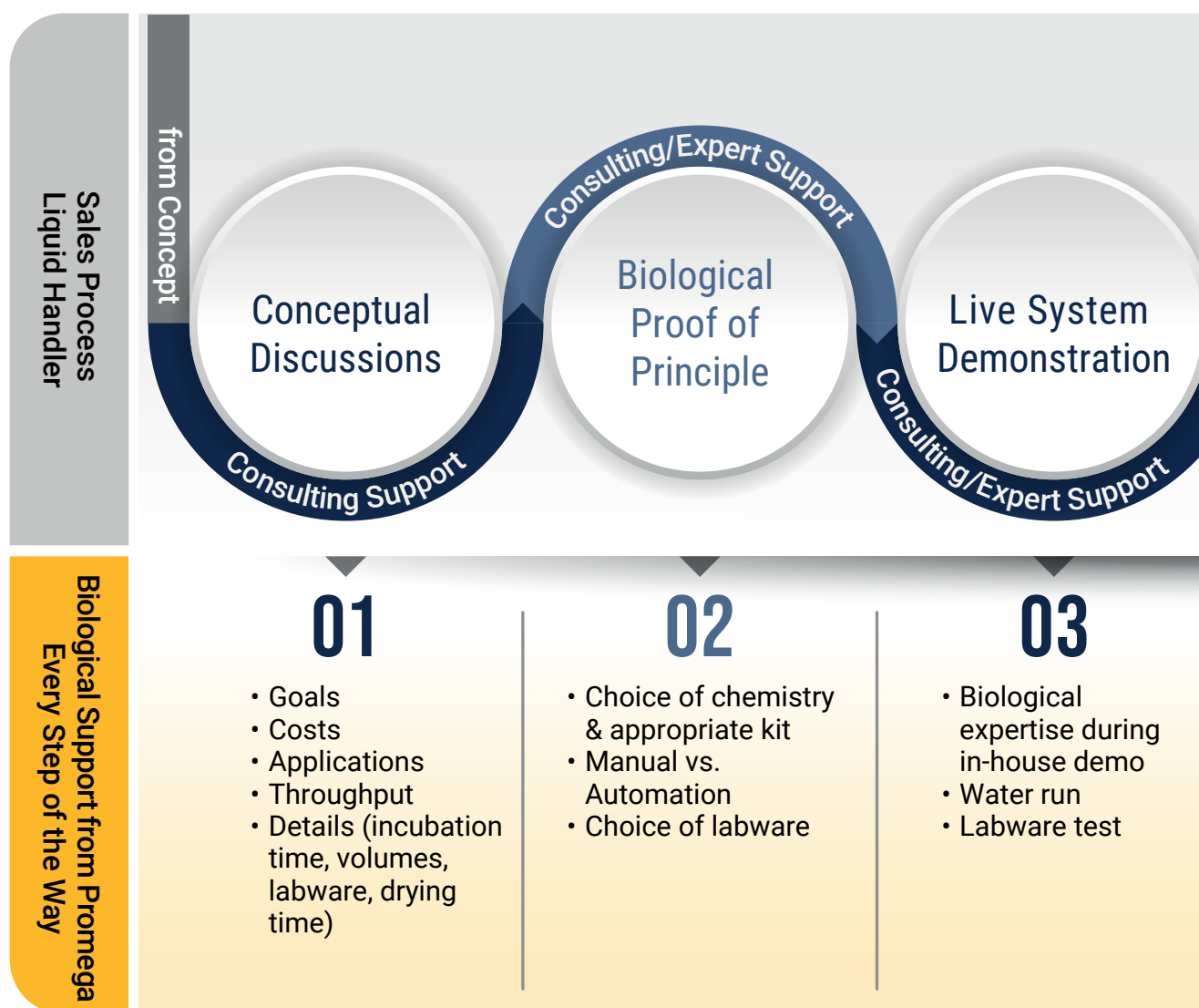
Cellular Analysis



- ▶ Cell Health Assays
- ▶ Cellular Metabolism Assays
- ▶ Immunoassays
- ▶ Bioassays

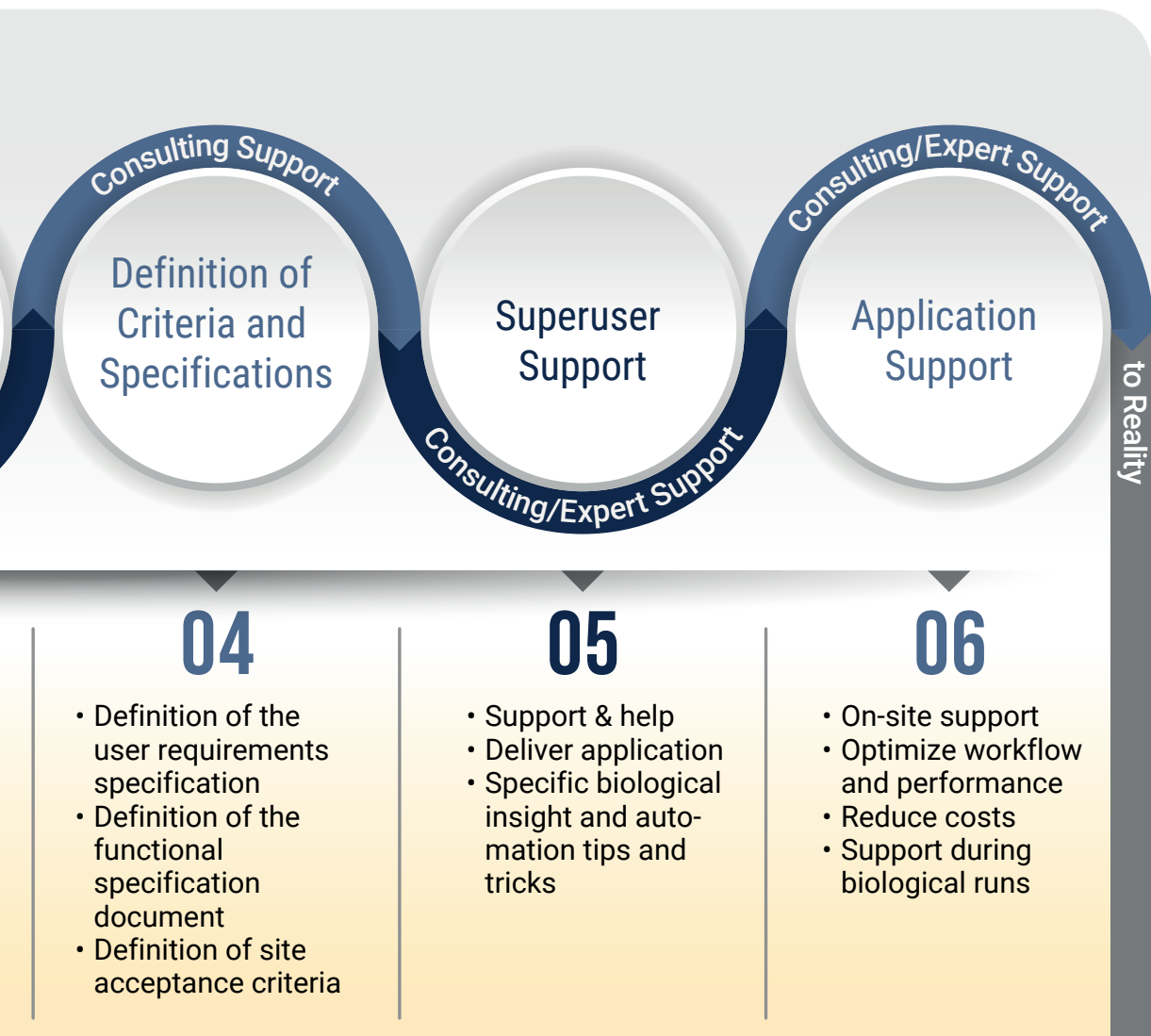
From Concept

The Automation Journey



From Concept to Reality and Beyond: Starting a lab automation project opens the door to more efficient, scalable, and reliable workflows. But moving from the first idea to a working solution can be complex. Success depends on a well-structured process: setting clear goals, understanding application needs, choosing the right tools and reagents, and aligning every element into a coordinated workflow.

to Reality

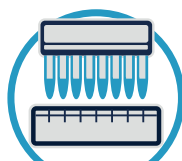


Throughout the journey, expert support helps ensure that each step—from initial discussions to system validation and biological applications—is carefully planned and executed. With Promega as a trusted partner, chemistry, instrumentation, and biology come together in a solution built around your specific needs. The result is a workflow that fits seamlessly into your lab, supported by a partner who remains by your side even beyond implementation.

Platforms



Ready-to-Go
Extraction Kits



Expert Support for Any
Liquid Handling System



Assay Scale-Up and
Custom Support

Experienced Partners

Streamlining your workflow becomes simpler with the dedicated support of an experienced partner. Our Field Support Scientists bring expertise across all major automation platforms, including:

ThermoFisher

KingFisher™ APEX &
KingFisher™ Flex

Analytik Jena

CyBio® FeliX

Hamilton

Microlab® STAR and STARlet

Eppendorf

epMotion®

Tecan

Fluent®, Veya™ and
Freedom Evo®

Beckman

Biomek® i5 and Biomek® i7

References

Plasmid Purification

High-Throughput Plasmid
Purification with Large-Volume
Bacterial Cultures on the
Hamilton STAR Liquid Handler



DNA Purification

Automated High-Throughput
DNA Purification from Blood
on the KingFisher™ Apex
PA929



RNA Purification

RNA Purification from Plant
Leaf and Seed using the
Maxwell® HT simplyRNA
Kit on the KingFisher™ Flex
Automation Platform PA130



RNA Purification

High Throughput Automated
RNA Purification From
Cultured Cells On The Tecan
Freedom EVO®150 PA648



Cell Viability Analysis

Automated CellTiter-Glo® 2.0
Cell Viability Assay on the
CyBio® FeliX from Analytik
Jena PA938



Compound Affect Testing

CyBio® FeliX flexible pipetting
platform with RealTime-
Glo™ MT Cell Viability Assay
multiplexed and CellTox™ Green
Cytotoxicity Assay PA933



Versatility

Selection of Sample Types for Nucleic Acid Purification



Selection of Products for Different Applications

Product Category	Product	Size	Cat. #
RNA Extraction from tissue	Maxwell® HT simplyRNA Kit, Custom	4 x 96 preps	AX7890
Genomic DNA Extraction	Maxwell® HT Genomic DNA Kit	4 x 96 preps	A6050
Plasmid Purification	Wizard MagneSil Tfx System	4 x 96 preps	A2380
ccfDNA Extraction	Maxwell® HT ccfDNA Kit	1 x 96 preps	A6030
Bioassay	PD-1/PD-L1 Blockade Bioassay	1 kit 2 x 1.0 ml	J1250 J1252
Cell Health	CellTiter-Glo® 2.0 Cell Viability Assay	10 ml 100 ml	G9241 G9242
RealTime-Glo™ MT Cell Viability Assay	RealTime-Glo™ MT Cell Viability Assay	100 reactions 1000 reactions	G9711 G9713

For further products & product information reach out to www.promega.com/support/tech-support

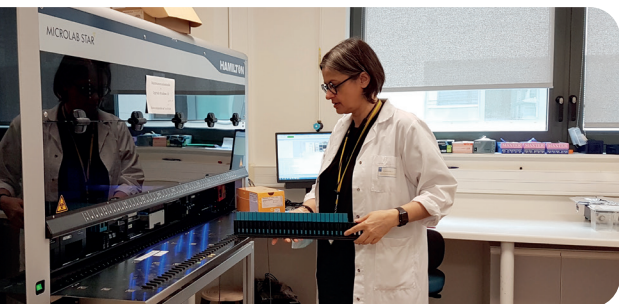
For more resources on lab automation, visit: www.promega.de/labautomation_resources

Legal Information: Trademarks: RealTime-Glo™ and CellTox™ are trademarks of Promega Corporation. KingFisher™ is a trademark of Thermo Fisher Scientific. Veya™ is a trademark of Tecan. Registered Trademarks: CellTiter-Glo® and Maxwell® are registered trademarks of Promega Corporation. Microlab® is a registered trademark of Hamilton Company. Fluent® and Freedom Evo® are registered trademarks of Tecan. Biomek® is a registered trademark of Beckman Coulter. epMotion® is a registered trademark of Eppendorf.

Customer

Streamlining Cancer Diagnostics

University Hospital of Rennes



“Promega team perfectly understood our needs and translated it into solutions. [...]. Nans [a Promega Field Support Scientist] came in November 2020 to install the script and test the protocol, between two [Covid] lockdown periods. The week after, I was already extracting FFPE samples with the new method. [...]. After we got the first results from samples extracted with the Promega Maxwell® HT FFPE System, we realized that sample quality with our in-house technique was suboptimal. Initially, ovarian tumor samples were tested with

Promega’s protocol then we extended its use to all tumor types. Quality improvement is quite phenomenal, we can reduce the number of re-testing and reduce time to results. [...]. We switched to routine on April 1st . [...]. After several weeks, we were able to make small adjustments to make everyone feel comfortable with the instrument and the protocol.”

Dr. Alexandra Lespagnol, Technical Manager

Background:

The Molecular Genetics of Cancer core-lab at the University Hospital of Rennes faced increasing demands for higher throughput and improved accuracy in nucleic acid extractions from challenging FFPE tissue samples. Lung cancer biopsies, often of low quality or limited in size, were especially problematic, as successful DNA extraction is critical for Next-Generation Sequencing (NGS) analysis. With their existing manual processes, the lab struggled to meet the standards required for reliable diagnostics and timely patient care.

Recognizing the need for automation, the lab partnered with Promega to develop a solution. The goal was clear: streamline the extraction process while improving both yield and quality of DNA from FFPE samples. Promega’s Field Support Scientist (FSS) team worked closely with the lab to implement the Maxwell® HT DNA FFPE chemistry on their existing Hamilton Star robotic platform. The result? A fully automated workflow capable of processing up to 96 samples in under two hours—vastly improving both efficiency and reproducibility.

Revolutionizing Plant Pathogen Detection

Valto Biocontrol

“The Maxwell® technology helped us isolate viral RNA of very high quality and high throughput. This made screening high numbers of samples easy and reliable. [...] . A huge benefit is that we can isolate both RNA and DNA or a mix using the same kit. [...] Promega’s Field Support Scientist team was open to converting protocols to be used on this isolation robot, from the custom dispensing of the consumables to a software file ready for the KingFisher itself.”

Dr. Menno Westerdijk, Head of Laboratories



Success Stories

Background:

Valto Biocontrol, an innovator in sustainable crop protection, needed a faster, more efficient way to detect plant pathogens like the Pepino mosaic virus. Their reliance on outsourced nucleic acid extraction resulted in delays of up to two weeks, creating bottlenecks that slowed decision-making and risked production setbacks for their V10 biocontrol product.

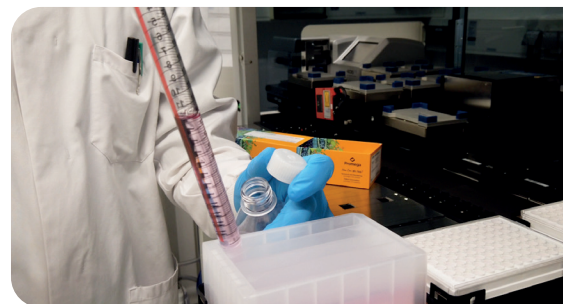
To take control of their workflow, Valto partnered with Promega to implement an automated extraction solution using the Maxwell® HT simplyRNA Kit. Promega's Field Support Scientist (FSS) team worked closely with Valto, developing a tailored high-throughput protocol to integrate seamlessly into their workflow. The automated solution allowed Valto to bring nucleic acid extraction in-house, cutting processing time from two weeks to just one day. It also improved data quality and reliability, strengthening their ability to manage pathogen detection and crop protection more effectively.

Automating Bioassays

Eurofins Biopharma Product Testing

"Working with Promega has been an excellent collaboration. Their in-depth knowledge of both their kits and the automation system itself has been invaluable. Our goal in this partnership is to implement a semi-automated and GMP-qualified cell-based potency assay (bioassay) that we can offer to our customers as a ready-to-use solution—without long delays or extensive method development. Promega's expertise and support are helping us bring this vision to life efficiently and seamlessly."

Dr. Sean Lin, Group Leader, Biologics and Bioassay Testing




Background:

Eurofins Biopharma Product Testing Munich GmbH (Eurofins Munich) faced an increasing demand for GMP sample testing to support regulatory-relevant potency assessments while ensuring method consistency, turnaround time and cost efficiency. However, their established workflows relied on manual method performance, which limited scalability and flexibility. To overcome these challenges, Eurofins Munich collaborated with Promega to implement an automated cell-based PD-1/PD-L1 Blockade Bioassay using the Hamilton Microlab STAR liquid handling system. A dedicated Promega Field Support Scientist worked on-site for two weeks with the Eurofins Munich Hamilton Specialist to program and implement the method.

This joint venture paved the way for GMP method qualification by increasing throughput and reducing hands-on time while maintaining method accuracy and precision comparable to manual methods. This advancement ensures compliance with industry regulations and expands Eurofins Munich's capacity to handle large sample volumes. Building on the success of this initiative, further semi-automation of other cell-based potency assays, such as ADCC and GLP-1 bioassays, is planned.

Contact

 Headquarters

 R&D and
Production

 Branches

 Distributors

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

Europe

Germany/Austria

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

Switzerland

Tel +41 44 878 90 00
ch_custserv@promega.com

France

Tel +33 0437 2250 00
Fax +33 0437 2250 10
fr.contact@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/Netherlands/Luxembourg

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

Spain/Portugal

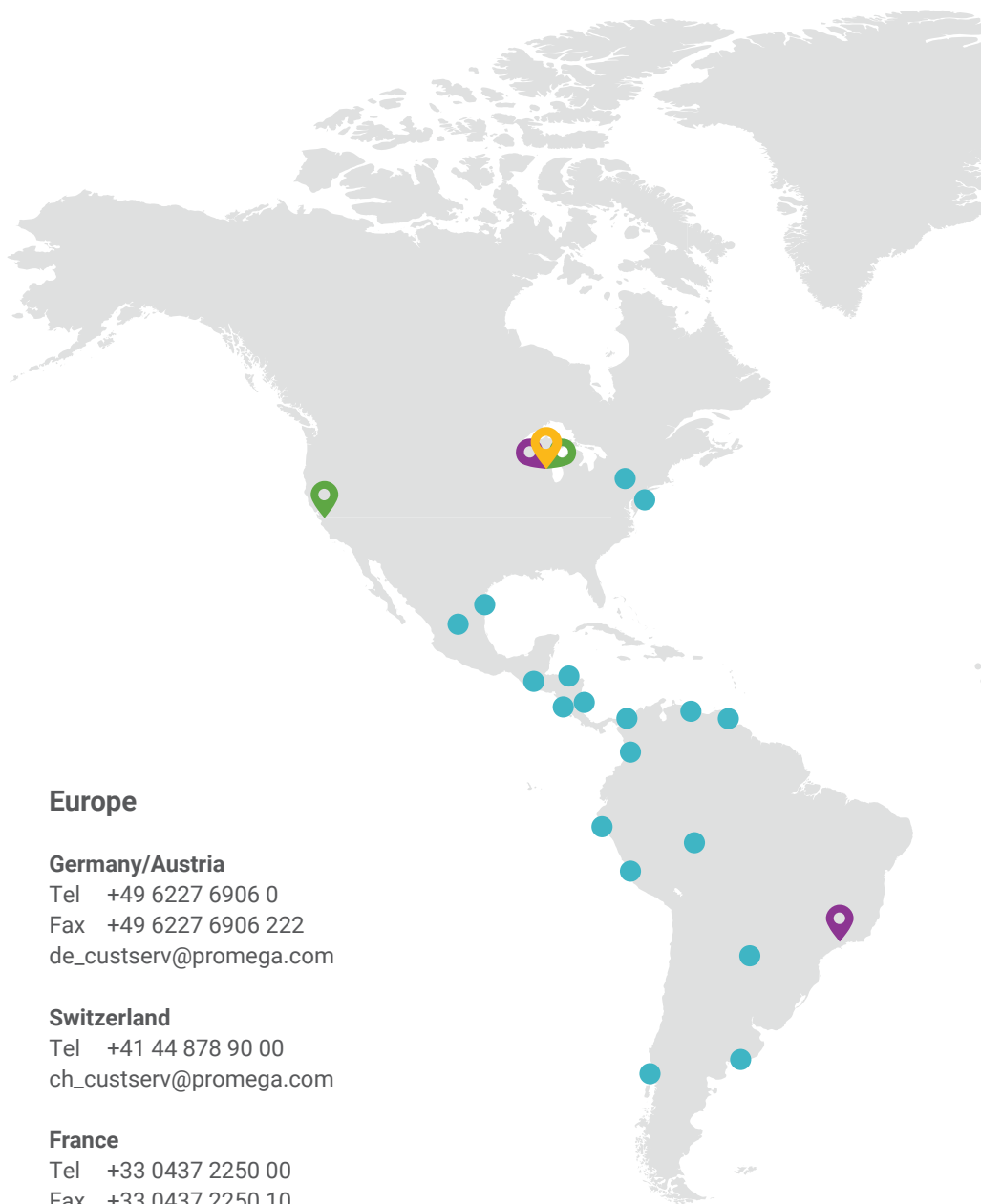
Tel +34 91 662 11 26
esp_custserv@promega.com

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





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

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 and Oceania

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)**

