



Promega

Promega Biotech AB

SE 020 402300 • sweorder@promega.com • promega.com/se
DK 800 8452 2450 • denorder@promega.com • promega.com/dk

Welcome to the Promega Life Science Catalogue 2008–2009. We are continually expanding our product line to meet your needs in a wide variety of scientific applications. In addition, we continue to update the format and indexing in our catalogue to help you find the products you need. Visit our website for online ordering, detailed product information and extensive support materials.

Ordering Made Easy

To place or enquire about your order please use one of the following convenient options:

PROMEGA BIOTECH AB

SWEDEN

Telephone: 020 402300

Customer Service Representatives are available Monday to Friday, 9:00 to 16:00

Online: promega.com/se/

Fax: 020 403880

E-mail: sweorder@promega.com

Mail: Promega Biotech AB
Finnboda Varvsväg 16
SE 131 73 Nacka

DENMARK

Telephone: 800 8452 2450

Customer Service Representatives are available Monday to Friday, 9:00 to 16:00

Online: promega.com/dk/

Fax: 800 8452 2455

E-mail: denorder@promega.com

Mail: Promega Biotech AB
Finnboda Varvsväg 16
SE 131 73 Nacka

To place your order, please be ready to provide the following information:

- Customer account number
- Name of researcher who will be using the product
- Purchase order number
- Catalogue number, product name and quantity of items you would like to order. In the event of a discrepancy the catalogue number will be used to fulfill the order
- Invoice address
- Shipping address, including department (if different)
- Your name and telephone number

Customized Products

—Buy exactly what you need

Promega offers custom services on all listed products targeting the areas of genomics, proteomics, cellular analysis and high throughput screening. Most standard catalogue products can be Customized to meet your specific application needs delivering exactly what you want for your research, improving your efficiency, saving you money and ensuring you get consistent reactions. Promega Custom offers:

- Novel formulations of existing products
- Custom concentrations
- Custom volumes
- Custom packaging and labeling
- Custom kit manufacturing
- Custom QC testing
- Manufacture of a “novel” product
- Custom protein purification

All Promega Custom products are designed and produced under the ISO 13485 quality standard and meet the same quality assurance standards as our catalogue products.

PromegaExpress™ Intelligent on-site product stocking programme

PromegaExpress™ provides the convenience of an on-site supply of freshly stocked Promega products to your lab giving you immediate access to a tailored range of products.

The programme offers the following advantages:

- Automatic, paper-free product replenishment and inventory management
- 24 hour, immediate access to a tailored range of Promega products using radio frequency (RF) identification, tagging and scanning
- Controlled access using unique pass key technology
- On-line inventory management system
- Product inventory customized to meet your requirements
- Durable, state-of-the-art freezers
- Dedicated support – Sweden 020 402300
Denmark 800 8452 2450

In addition, we still run the standard PromegaExpress™ programme offering on-site stock stored in standard lockable freezers or cupboards for room temperature storage. Both systems are backed by comprehensive support from Promega to ensure the products meet your research requirements.

For further information about installing PromegaExpress™ please contact our dedicated PromegaExpress™ support team in Sweden 020 402300 or in Denmark 800 8452 2450. On-site freezer installation is subject to meeting certain business conditions.

Pricing/Terms

Prices are subject to change without notice. Payment terms are 30 days from invoice date.

Note: “Pls. Enq.” listed as the price means “please ask” about the price and/or availability of the product.

Bulk Orders

We are pleased to offer unpublished higher discounts on large volume purchases. Please contact the customer services team.

Payment Options

Please remit your payment via:

Org nr/VAT No SE556717789301

Bankgiro: 211-0781

Plusgiro: 31720-6

Bank: Nordea, 105 71 Stockholm

Account No: 3201 17 10979

BIC/SWIFT: NDEASESS

IBAN: SE873000000032011710979

Confirming Orders

To make order processing quicker and more efficient, Promega does not require a written confirmation of product orders. If you choose to place an order in writing, we ask that you do not telephone us with the same order, since this would result in a duplicate order.

To avoid duplication, please ensure that any written confirmation of a telephone, fax, e-mail or online order is clearly marked “CONFIRMATION”.

Delivery

Orders received prior to 16:00 Monday to Wednesday are delivered within 48 hours (standard delivery).

Orders received on Wednesday are delivered on Friday, subject to Institution receiving hours.

Orders with a value less than 3500 SEK are subject to 150 SEK standard delivery charge. Orders with a value greater than 3500 SEK have free of charge standard delivery (48 hours).

We offer 24 hour delivery, subject to premium surcharge of 150 SEK. Dry ice deliveries are subject to 80 SEK surcharge

If you need to know when your order has been delivered, we can provide notification of delivery on request. Simply provide us with your email address and ask for a delivery notification when you place your order.

Returns

Because of the high sensitivity of our products we are not able to accept all returns. Our commitment to product quality and performance prevents us from reselling returned goods. If you would like to initiate a return, please contact Promega to obtain a return authorisation code. In instances of purchasing errors, we are unable to accept the return of products.

Recyclable Packaging

Promega strives to use packaging material that maximizes product integrity and performance, while minimizing impact on the environment. All products incorporate highly recyclable packaging materials. Product systems are packaged in paper-based, non chlorine bleached cardboard, coated for convenient opening when stored in the freezer.

We ship orders of heat labile products in EPS (Expanded Polystyrene) foam boxes. We believe that EPS foam is the best available product for temperature maintenance, cushioning and durability.

The Promega Points Reward Scheme

Every time you order a product from Promega we automatically include Promega Points for you to collect to redeem for one of the great rewards we have on offer including:

- Gift vouchers
- Lab equipment
- Journal subscriptions
- Thesis binding support
- Books
- Fun Stuff

For further information go to:

promega.com/se/swecollection.htm

Electronic Procurement

Promega is dedicated to improving your procurement processes and reducing transaction costs.

Our eBusiness capabilities extend from Procurement through Payment including:

- Hosted catalogs
- PunchOuts
- XML ordering/acknowledgement/notification
- Electronic invoicing

Please contact our eBusiness Team at b2b@promega.com to discuss your requirements.

Promega Biotech AB has e-invoicing capabilities. Please contact customer services.

Technical Services

Promega has Ph.D. qualified Technical Services Scientists equipped with the lab experience vital for responding to your enquires. You are our major source of information about product performance, new lab applications and helpful lab hints. We actively encourage you to contact us using any of the following convenient methods:

Technical Services Scientists are available Monday to Friday from 9:00 to 16:00

SWEDEN

Telephone: 020 402300

Online: promega.com/techserv/

Fax: 020 403880

E-mail: swetech@promega.com

Mail: Promega Biotech AB
Finnboda Varvsväg 16
SE131 73 Nacka

DENMARK

Telephone: 800 8452 2450

Online: promega.com/techserv/

Fax: 800 8452 2455

E-mail: dentech@promega.com

Mail: Promega Biotech AB
Finnboda Varvsväg 16
SE131 73 Nacka

Technical Sales Specialists (TSS)

Promega's Technical Sales Specialists (TSS) serve as the principal representatives of the Promega technical sales programme in a specified geographic territory. Each TSS has solid experience and training in the processes and techniques employed in today's research laboratories.

Promega On Line

promega.com

In addition to global news and announcements, the Promega web site is continually pushing ahead to provide you with improved internet features, like our online ordering system and the best technical product support in the industry.

For information specific to Promega Sweden: promega.com/se/ Denmark: promega.com/dk/

For everything technical about our products, look on the Promega home page for "Technical Resources" or go there directly at the following address: promega.com/techserv/

On the Technical Resources page, you will find links to a range of information including:

- Full text & graphic versions of all: Technical Bulletins and Manuals (promega.com/tbs/); and, Vector Sequences (promega.com/vectors/).
- Complete articles from current: eNotes (promega.com/enotes/); Promega Notes (promega.com/pnotes/); Cell Notes (promega.com/cnotes/); and, Profiles in DNA (promega.com/profiles/) magazines.
- Technical tools such as: Citations, calculators, transfection assistants, tools for cell-based assays, etc.
- Special Offers: Visit Promega AB for exceptional offers (promega.com/se/offers)



TERMS AND CONDITIONS OF SALE

BASIS OF SALE

Promega Biotech AB ("Promega") will sell and the person, firm, company or institution ("Buyer") will purchase the goods listed on Promega's current product and price list ("Goods") subject to these terms and conditions.

No quotation given by Promega shall constitute an offer to supply the Goods. An order placed by the Buyer, whether or not in response to a quotation, shall constitute an offer made to Promega subject to these terms and conditions. No terms or conditions put forward or implied by the Buyer, in the order, correspondence or elsewhere shall be binding on Promega, such conditions or stipulations are hereby excluded and extinguished.

All Contracts for Goods are subject to availability.

Any error or omission in any quotation, catalogue, sales literature, invoice or other document issued by Promega shall be subject to correction by Promega without liability to Promega.

ORDERS

The buyer shall be responsible to Promega for the accuracy of all orders. In the event of a discrepancy between catalogue number and its product description, the catalogue number will be used to satisfy the order.

The buyer shall be responsible to Promega for all confirmation orders to be clearly marked CONFIRMATION to avoid duplication of an order.

All Buyer's orders are subject to Promega's acceptance.

Promega reserves the right to make changes in the specification of Goods ordered which do not materially affect quality or performance of the Goods.

PRICE

Unless specifically agreed otherwise, the price for the Goods will be that as shown in Promega's current catalogue at the time of delivery.

Promega reserves the right without notice to vary or amend the price for the Goods shown in any price list, literature or in any other document issued by Promega. Such price changes shall not apply to accepted orders.

The price is exclusive of Value Added Tax and all other similar sales taxes.

Promega reserves the right to charge a delivery, packaging and/or handling charge on all orders.

PAYMENT

Our terms are strictly 30 days from date of invoice. No settlement discounts or other deductions may be made against amounts due on invoices.

The time for payment shall be of the essence for all contracts for Goods. In the event of default in payment by a due date, Promega reserves the right, without prejudice to other rights, to: (i) charge interest at the base rate +2% as of the time of default and/or (ii) suspend further deliveries.

For Buyers with an agreed credit account facility with Promega, unless specifically agreed otherwise, the payment terms are 30 days from invoice date.

For Buyers without an agreed credit account facility with Promega, unless specifically agreed otherwise, Promega is to have received cleared funds for all amounts due before dispatch of the Goods.

All amounts due are payable in Swedish Kronor or any lawfully recognized currency that may replace the Swedish Kronor by draft on a Swedish bank unless otherwise agreed in writing.

All costs including legal expenses incurred by Promega in recovering overdue debts will be payable by the Buyer.

Promega reserve the right to withhold further supplies in the event of amounts payable being overdue, breach of any of these conditions or for any other reason which at the discretion of Promega warrants such action.

DELIVERY AND RETURN

Delivery of Goods shall be at Buyer's cost, to the Buyer's place of business as notified in writing to Promega. Delivery of Goods shall be at the Buyer's place of business as notified in writing to Promega. Applicable delivery charges are pre-paid by Promega and added to the Buyer's invoice as a separate line item.

Goods supplied will not be accepted for return unless returned in accordance with an authorisation issued by Promega.

RISK AND PROPERTY

Risk shall pass to the buyer:

Where Goods are stored by Promega upon such Goods being stored; or in all other cases upon delivery or attempted delivery (where Buyer refuses delivery).

Property in the Goods shall not pass to the Buyer until all amounts due have been received by Promega in full.

WARRANTIES AND LIABILITIES

Subject as mentioned below. Promega warrants that its Goods perform as described in the Promega product information sheets, which state the conditions and results of product performance tests conducted in our quality assurance laboratory. These results are warranted to meet or exceed our published specifications when used under normal conditions in a laboratory.

To the fullest extent permitted by law, Promega's total liability to the Buyer under these terms and conditions is subject to the conditions of this Section.

Should any product fail to perform as warranted, Promega's sole liability shall be limited to such replacement of the Goods or full credit of the purchase price at the discretion of Promega.

The above warranty is subject to:

Promega being under no liability for any defect in the Goods resulting from Buyers designs or requirements, willful damage, negligence, failure to follow instruction, misuse, use beyond its expiry date (if applicable) or if the amounts due have not been paid in full to Promega.

Promega being under no liability for any incidental, consequential or contingent damages.

All other warranties and conditions expressed or implied are expressly excluded to the fullest extent permitted by law.

HAZARDS AND TOXICITY

All Goods offered by Promega are intended for use by qualified professionals who are trained in good laboratory procedures and are familiar with potential hazards.

Promega Goods having known toxicity are shipped with a Material Safety Data Sheet which describes, to our knowledge, the potential dangers. The absence of a toxicity warning does not, however, preclude a possible health hazard.

PRODUCT USE LIMITATIONS, TRADEMARKS, SERVICE MARKS AND PATENTS

Unless specified otherwise, all Promega goods are sold for research and/or in vitro use only.

Please see Technical and Legal References section of this current catalogue for acknowledgements, disclaimers, trademarks, service marks and patent information.

GENERAL

Headings are for reference only and shall not affect interpretation.

Notices shall be served by post on the registered offices of the parties or principal place of business.

No waiver of a breach of contract by Promega shall operate as a waiver of any other or subsequent breach. If any provision of these conditions is held invalid or unenforceable in whole or part, the validity or enforceability of the other provisions shall not be affected.

These Terms and Conditions and any Contract are governed by Swedish law. Any dispute arising out of or in connection with these Terms and Conditions shall be determined by the Swedish courts.

QUALITY MANAGEMENT SYSTEM COMMITMENT— ISO 13485

Promega Corporation has a long history in supporting its customers with high-quality products, services and information. Promega first certified to international standards for quality management systems in 1998, and this commitment continues with our current ISO 13485-2003 certification, required for the development, manufacture, testing and delivery of medical devices around the world. The ISO series of quality management system standards are developed and maintained by the International Organization for Standardization. An organization achieving ISO certification has demonstrated to a third party body of experts that the organization meets all requirements of the standard. This certification assures our global customers that Promega has established reliable and effective processes for product development, manufacturing, distribution and customer support in its quality management system. With this certification, we establish our commitment to our customers, to our business, and to all those who rely on and benefit from the use of our products.

Customers may obtain a copy of Promega's current ISO 13485 Certificate at: promega.com/iso/

