

Terms and Conditions of Sales

Ordering

You have the following options when ordering products directly from Jena Bioscience:

- Mail orders
- Telephone orders
- 24 hour fax ordering
- Online ordering

Please provide us with the following information when ordering:

- Your name, name of institution
- Billing and shipping address
- PO number (if applicable)
- Catalog number of products and quantity needed
- Contact person and contact data for questions

Mail orders

Please send your mail orders to the following address:

Jena Bioscience GmbH
Loebstedter Strasse 80
07749 Jena
Germany

Telephone orders

We will accept telephone orders from Monday to Friday between 8:00 am and 16:00 pm Central European Time.
+49 – 3641 – 628-5000

24 hour fax ordering

Please send your fax order to:
+49 – 3641 – 628-5100

Online ordering

Jena Bioscience products can be ordered online. When ordering by e-mail, please direct your orders to
orders@jenabioscience.com

Products can also be ordered online in our online shop:
<http://www.jenabioscience.com>

Important Notice:

Products that have been ordered by mistake cannot be returned to Jena Bioscience. Products that are returned unrequestedly to Jena Bioscience will not be accepted, but fully charged to the customer's account.

Jena Bioscience GmbH

Loebstedter Str. 80
07749 Jena, Germany
Phone: +49 – 3641 – 628-5000
Fax: +49 – 3641 – 628-5100
Email: info@jenabioscience.com
<http://www.jenabioscience.com>

Management:

Thomas Billert, COO
Dr. Mathias Gruen, CEO
VAT no.:DE 195825742
Amtsgericht Jena
HRB 207171

Bank Data:

Bayerische Hypo- und Vereinsbank AG Jena
Account no.: 419 60 90
Bank Code : 830 200 87
IBAN: DE 05830200870004196090
SWIFT: HYVEDEMM463

Shipping

All customers will receive a fax confirmation of the order with invoice and shipping waybill number.

International orders are shipped either by General Overnight, by FedEx or by UPS Express service, depending on the customer's location and on the products to be shipped. Domestic shipments within Germany are sent by General Overnight Express service. If you wish your order to be shipped by a different carrier, please contact us and provide all necessary information with your order.

All orders are shipped EXW (Incoterms 2000). Please contact us if a different shipping term is required for your order.

Prices and Charges

Please note that the prices of products in the catalog and on our website do not include freight charges, duties, taxes or customs fees.

Freight charges will be prepaid and added to the invoice. Freight charges for online orders are indicated when you check out of the online store. If you need information on freight charges for your particular order, please contact us with all necessary information.

Jena Bioscience will not pay any duties, taxes or customs fees.

Products and prices are subject to change without notice. Current pricing will be confirmed at the time of your order. No minimum order required.

We offer free shipping of all orders worth EUR 250+ (excl. VAT) within Germany.

Payment

Invoices will be issued after your order has been shipped and will be sent to the billing address by separate mail. Invoices will not be included within the shipments. In case of partial deliveries, separate invoices will be issued after each shipment has left Jena Bioscience. You will find payment information (bank addresses and account data) on each invoice. Jena Bioscience accepts payment by:

Check

Please send your payment checks to the following address:

Jena Bioscience GmbH
Loebstedter Strasse 80
07749 Jena
Germany

We kindly ask you to make sure that our invoice number and your customer number appear on the cheque.

Jena Bioscience GmbH

Loebstedter Str. 80
07749 Jena, Germany
Phone: +49 – 3641 – 628-5000
Fax: +49 – 3641 – 628-5100
Email: info@jenabioscience.com
<http://www.jenabioscience.com>

Management:

Thomas Billert, COO
Dr. Mathias Gruen, CEO
VAT no.:DE 195825742
Amtsgericht Jena
HRB 207171

Bank Data:

Bayerische Hypo- und Vereinsbank AG Jena
Account no.: 419 60 90
Bank Code : 830 200 87
IBAN: DE 05830200870004196090
SWIFT: HYVEDEMM463

Wire transfer

Please remit your payments to one of the following bank accounts:

HypoVereinsbank - UniCredit Bank AG	Account No.: 4196090
Niederlassung Thüringen	Bank code (BLZ): 83020087
Schillerstrasse 4	IBAN: DE
07745 Jena	05830200870004196090
Germany	SWIFT: HYVEDEMM463

Sparkasse Jena-Saale-Holzland	Account No.: 32417
Ludwig-Weimar-Gasse 5	Bank code (BLZ): 83053030
07743 Jena	IBAN: DE 22830530300000032417
Germany	SWIFT: HELADEF1JEN

Credit card

Jena Bioscience accepts the following credit cards:

- VISA
- Mastercard
- American Express



If you wish to pay by credit card, please provide the following credit card information:

- Card number
- Expiry date
- Security code (VISA/Mastercard: 3 digits, to be found on your card's back side in the upper right corner of the signature field; AmEx: usually 4 digits (sometimes only three), to be found on the front side of your card above the card number)

Patent Disclaimer

Unless explicitly stated, no license or immunity under any patent is either granted or implied by the sale of any of our products. Jena Bioscience does not warrant that the resale or use of its products delivered will not infringe the claims of any patent, trademark or copyright covering the use of the product itself or its use in the operation of any process. Furthermore, the purchaser assumes all risks of patent, trademark or copyright infringement associated with any such use, combination or operation.

Jena Bioscience GmbH

Loebstedter Str. 80

07749 Jena, Germany

Phone: +49 – 3641 – 628-5000

Fax: +49 – 3641 – 628-5100

Email: info@jenabioscience.com

<http://www.jenabioscience.com>

Management:

Thomas Billert, COO

Dr. Mathias Gruen, CEO

VAT no.:DE 195825742

Amtsgericht Jena

HRB 207171

Bank Data:

Bayerische Hypo- und Vereinsbank AG Jena

Account no.: 419 60 90

Bank Code : 830 200 87

IBAN: DE 05830200870004196090

SWIFT: HYVEDEMM463