

**Please fill in your order and fax it to:
+49 – 3641 – 628 5100
Jena Bioscience Fax Order Form**

Billing address:

Name

University/Company

Institute/Department

Address

Address

Postcode, City, Country

Telephone

Fax

Email

Shipping address:

Name

University/Company

Institute/Department

Address

Address

Postcode, City, Country

Telephone

Fax

Email

Customer number

Your PO number

VAT number (EEC customers only)

Date, Signature

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



Card holder

Expiry date

Card number

Security code
(VISA/Mastercard: 3 digits on card's back side, upper right corner of signature field;
AmEx: 4 digits, card's front side, above card number)

	Catalog number	Product	Quantity	Net Price per Item EURO	Net Price all Items EURO
1					
2					
3					
4					
5					
6					
7					
8					
9					
10					
				Total	

Jena Bioscience GmbH
Loebstedter Strasse 80
07749 Jena, Germany

Tel.: +49 – 3641 – 628 5000
Fax: +49 – 3641 – 628 5100
<http://www.jenabioscience.com>