

Please copy this page, fill in your order and fax it to:
+49 – 3641 – 628 5100

Jena Bioscience Fax Order Form

Shipping address:

Billing address:

Name

University/Company

Institute/Department

Address

P.O.Box

Postcode/City

Phone

Fax

Email

Customer number

Name

Address

Address

Address

Address

VAT number (EEC 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					
11					
12					
13					
14					
				Total	

Jena Bioscience GmbH
 Loebstedter Strasse 80
 07749 Jena, Germany

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