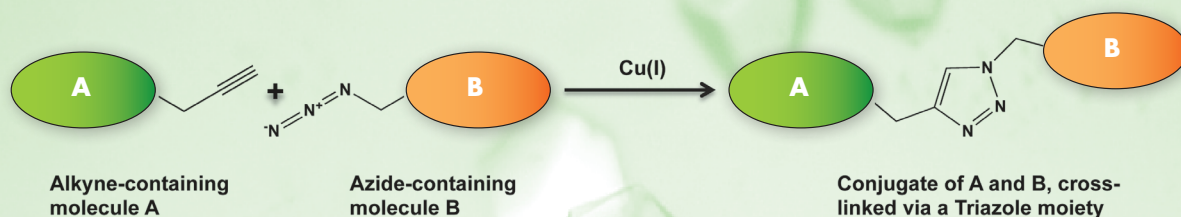


**NEW**

Click Chemistry Reagents

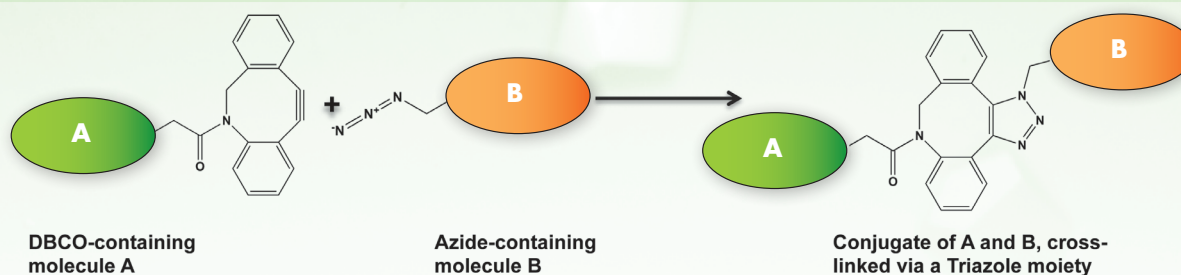
Click Chemistry has evolved into a state-of-the-art method for labeling and crosslinking of biomolecules. It combines selectivity, quantitative yields, and bio-orthogonality with working under physiological conditions. **Jena Bioscience offers a broad variety of reagents for click chemistry, including:**

Alkyne-Reagents for Cu(I)-catalyzed Click Reactions



- Click Chemistry Kits
- Alkyne-containing Nucleosides
- Alkyne-containing Nucleotides
- Alkyne-containing Phosphoramidites
- Alkyne-containing CPGs
- Alkyne-containing Spacers
- Alkyne-containing Fluorescent Dyes
- Auxiliary Reagents

DBCO-Reagents for Copper-free Click Reactions



- DBCO-containing Chemical Modification Reagents
- DBCO-containing Biotinylation Reagents
- DBCO-containing Spacers
- DBCO-containing Fluorescent Dyes

Azide-Reagents for Click Reactions

- Non-Fluorescent Azides
- Biotin Azides
- PEG and Spacer Azides
- Azides of Amino Acids
- Azides of Fluorescent Dyes
- Azides of Quencher Dyes
- Nucleotide Azides

Check out our complete click chemistry product portfolio and find more infos at

www.jenabioscience.com/click



Jena Bioscience GmbH
Loebstedter Str. 80
07749 Jena
Germany

Phone +49 (0)3641-62 85 000
Fax +49 (0)3641-62 85 100
www.jenabioscience.com
info@jenabioscience.com

 Jena Bioscience

Jena Bioscience Fax Order Form

Please copy this page, fill in your order and fax it to: **+49(0)3641-62 85 100**

Shipping address

Billing address

Name	Customer number
University / Company	University / Company
Institute / Department	Institute / Department
Address	Address
Postcode	Postcode
City / Country	City / Country
Phone	VAT number (EEC only)
Fax	PO number
E-mail	Date / Signature

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



Card holder Card number

Expiry date 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
1			
2			
3			
4			
5			
6			
7			
8			
9			
10			
11			
12			
13			
14			
15			



Jena Bioscience GmbH
Loebstedter Str. 80
07749 Jena
Germany

Phone +49 (0)3641-62 85 000
Fax +49 (0)3641-62 85 100
www.jenabioscience.com
info@jenabioscience.com

