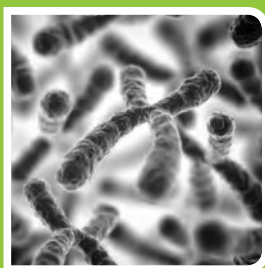


Probes: Non-radioactive Labeling



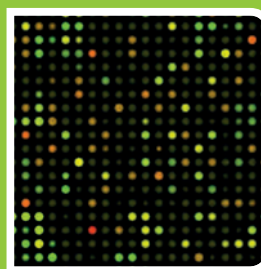
Historically, mainly radioisotope labels were used for biomolecule (e.g. DNA, RNA, protein) labeling however, in recent years **non-radioactive biomolecule labeling has become an attractive alternative over radioactive approaches**. This is due to significant improvements of the available detection methods (speed, sensitivity, and versatility) that make the **analytical performance of non-radioactive labels e.g. fluorescent dyes or haptens^[1] comparable to that of their radioactive counterparts**. In addition, **non-radioactively labeled biomolecules are clearly superior in many practical aspects** such as enhanced stability, convenient handling and greatly improved safety profiles.

DNA/cDNA Labeling



Fluorescent, Hapten^[1]-, Click Chemistry- and Amine- modified Nucleotides & Labeling Kits for the enzymatic preparation of labeled DNA/cDNA probes

RNA/cRNA Labeling



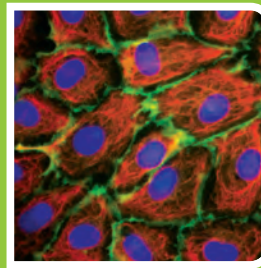
Fluorescent, Hapten^[1]-, Click Chemistry- and Amine-modified Nucleotides & Labeling Kits for the enzymatic preparation of labeled RNA/cRNA probes

Protein Labeling



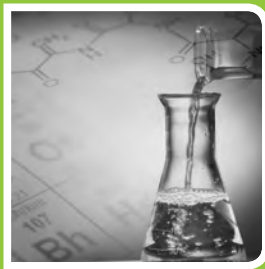
Novel & Classic Labeling Substances and Kits for Fluorescent, Hapten^[1]- and Luminescent-labeling of various protein sources, both random and site-directed (N- and/or C-terminal)

Cell Labeling



Fluorescent reporter and Protein & Nucleic Acid Internalization Kits

Reactive Components



Reactive Fluorescent, Hapten^[1]-modified, Luminescent and bifunctional reagents for labeling and crosslinking of biomolecules such as DNA, RNA and proteins

More Information



www.jenabioscience.com/probes

^[1] Haptens = Biotin, Digoxigenin, Dinitrophenol

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			