



#### About us

Jena Bioscience GmbH was founded in 1998 by a team of scientists from the Max-Planck-Institute for Molecular Physiology in Dortmund. 25+ years of academic knowhow were condensed into the company in order to develop innovative reagents and technologies for the life science market. Since the start up, the company has evolved into an established global reagent supplier with about 8000 products in stock and a customer base in 80+ countries. Jena Bioscience serves three major client groups:

- Research laboratories at universities, industry, government, hospitals and medical schools
- Pharmaceutical industry in the process from lead discovery through pre-clinical stages
- Laboratory & diagnostic reagent kit producers and re-sellers

Our company premises are located in the city of Jena / Germany.



Our company premises are located in the Saalepark Industrial Estate in the northern part of the city of Jena / Thüringen / Germany. In March 2015 we moved all operations to our own, new 2.500 som company building.



Jena Bioscience's products include nucleosides, nucleotides and their nonnatural analogs, recombinant proteins & protein production systems, reagents for Click Chemistry, for crystallization of biological macromolecules and tailormade solutions for molecular biology and biochemistry.

In our chemistry division, we have hundreds of natural and modified nucleotides in stock. In addition, with our pre-made building blocks and in-house expertise we manufacture even the most exotic nucleotide analog from mg to kg scale.

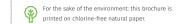
In the field of recombinant protein production, Jena Bioscience has developed its proprietary LEXSY technology. LEXSY (Leishmania Expression System) is based on a S1-classified unicellular organism that combines easy handling with a eukaryotic protein folding and modification machinery including mammalian-like glycosylation. LEXSY is primarily used for the expression of proteins that are expressed at low yields or are inactive in the established systems, and expression levels of up to 500 mg/L of culture were achieved.

For the crystallization of biological macromolecules – which is the bottleneck in determining the 3D-structure of most proteins – we offer specialized reagents for protein stabilization, crystal screening, crystal optimization and phasing that can reduce the time for obtaining a high resolution protein structure from several years to a few days.

Our reagents are complemented with a large selection of molecular biology kits for PCR, mutagenesis, cloning and alike as well with innovative reagents for the functionalization and labeling (fluorophores, haptens) of biomolecules with focus on Click Chemistry approaches.

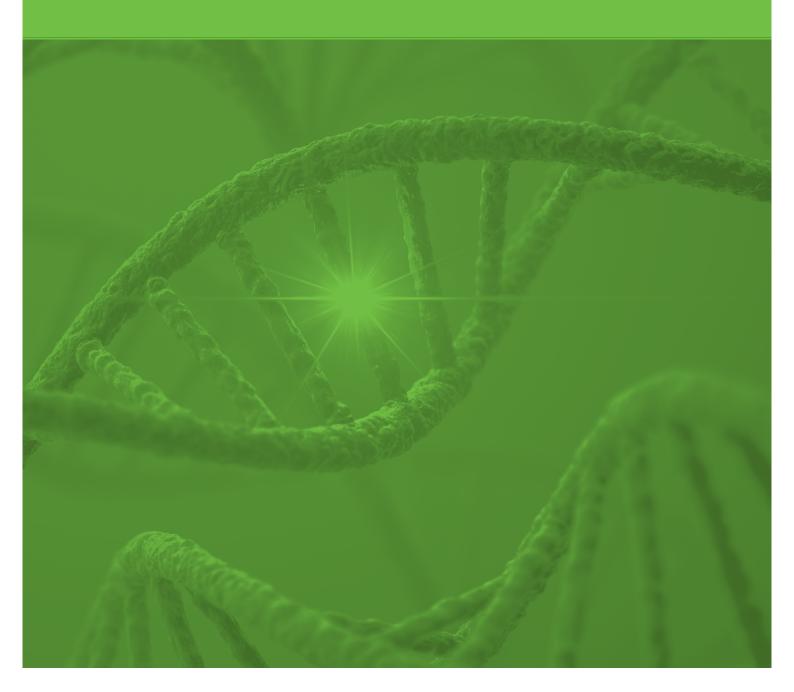
We combine highest quality standards for all our products (certified according to DIN EN ISO 9001) with individualized customer support. We establish direct lines of communication from clients to our in-house scientists, resulting in productive interactions among people with similar research interests who speak the same language. Furthermore, we offer support programs and attractive discount schemes for young scientists establishing their own labs. If you wish to receive more information, just send an e-mail to info@jenabioscience.com.





# **Table of contents**

Introduction	4
Labels and their detection	5
Incorporation of labels into DNA/RNA	6







## Introduction

DNA and RNA fragments of variable length ("Probes") are able to form non-covalent, highly specific duplexes with a complementary nucleic acid strand (termed hybridization). When a label is attached to such a hybridization probe it can thus efficiently serve for detection of a defined DNA or RNA target sequence.

Typical applications of hybridization probes are (fluorescence) in situ hybridization ((F)ISH), microarray-based gene expression profiling or electrophoretic mobility shift assays (EMSA) that are routinely used in research or clinical diagnostics to

- detect and localize specific DNA sequences and their potential aberrations (mutations, deletions, duplications) on chromosomes ((F)ISH, Fig. 1A).
- identify and analyze protein-DNA/-RNA interactions (EMSA, Fig. 1B).
- simultaneously measure the relative abundance of multiple RNAs (e.g. microarray-based gene expression profiling).

Historically, mainly radioisotope labels were used for hybridization probes however, in recent years the incorporation of non-radioactive labels has gained increasing popularity. This is due to significant improvements of the available detection methods (speed, sensitivity, and versatility) that make the analytical performance of non-radioactive labels comparable to that of their radioactive counterparts. In addition, non-radioactive probes are clearly superior in many practical aspects such as **enhanced stability, convenient handling and greatly improved safety profiles**.

Therefore, and together with the corresponding detection systems and well-established incorporation methods, non-radioactive nucleic acid labeling probes have become a superior alternative over radioactive approaches.

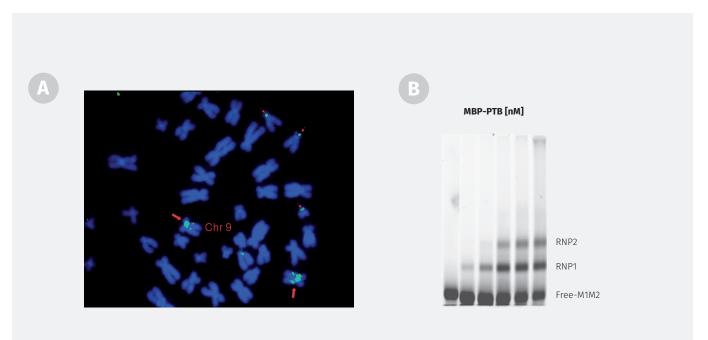


Figure 1

A: Chromosome 9 was visualized by fluorescence in situ hybridization (FISH). A DNA probe (Bac268E1) was fluorescently labeled by nick translation using Jena Bioscience's Aminoallyl-dUTP-XX-ATTO-488 (Cat.-No. NU-803-XX-488).

B: Polypyrimidine tract-binding protein (MBP-PTB) as a new component of the oskar mRNA/ ribonucleoprotein (M1M2/RNP1/2) complex was detected by an electrophoretic mobility shift assay (EMSA). The mRNA probe (M1M2) had been labeled by T7 RNA polymerase mediated in vitro transcription using Jena Bioscience's Aminoallyl-UTP-ATTO-680 (Cat.-No. NU-821-680) (adapted from [1]).

### Labels and their detection

The most commonly used labels for the generation of non-radioactively DNA or RNA hybridization probes are fluorophores and haptens, the latter meaning Biotin and Digoxigenin.

Fluorescent probes are detected directly after incorporation by fluorescence spectroscopy (Fig. 2A). Nowadays, a plethora of fluorescent dyes is available with optical properties that cover the whole UV-Vis spectrum and fit to common commercial light sources and filter systems. The dyes differ, however, greatly in terms of incorporation efficiency, photostability, hydrophilicity, and size/bulkiness. Classical dyes such as rhodamine, fluorescein and cyanine derivatives have been the most widely used so far but recently an increasing number of new dyes (e.g. ATTO dyes) has become very popular.

In contrast to fluorophores, Biotin and Digoxigenin are indirect labels since their visualization requires a secondary reporter molecule (Fig. 2B). The detection of Biotin, also known as vitamin H,

relies on its high affinity (KD= 10-15 M) to Streptavidin, a 60 kDa tetrameric protein purified from the bacterium *Streptomyces avidinii* [2]. Biotinylated hybridization probes are therefore easily detected through Streptavidin carrying either a reporter enzyme or a fluorescent dye (Fig. 2B.1). Since however, endogenous biotin from biological samples often interferes with biotin detection (resulting in high backgrounds and false positives [3][4]), the use of Digoxigenin, a steroid exclusively present in Digitalis plants, may sometimes be preferable [3][5]. Visualization of incorporated Digoxigenin is achieved in a similar manner by Digoxigenin antibody conjugates with a reporter enzyme or a fluorescent dye (Fig. 2B.2).

The choice of label and method of detection is generally determined by the particular downstream application. While speed of detection and ease of quantification are major advantages of fluorescent labels, indirect methods with secondary reporter molecules allow signal amplification with resultant increased sensitivity.

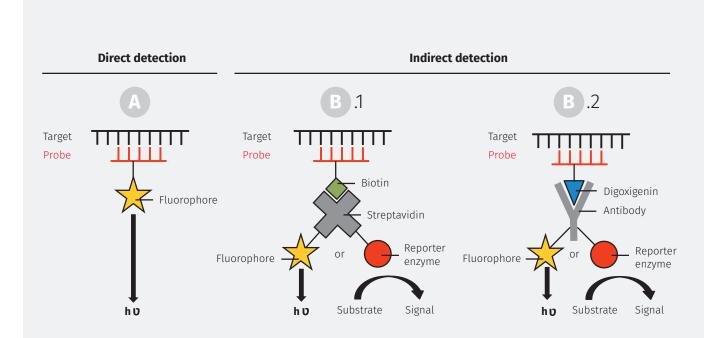


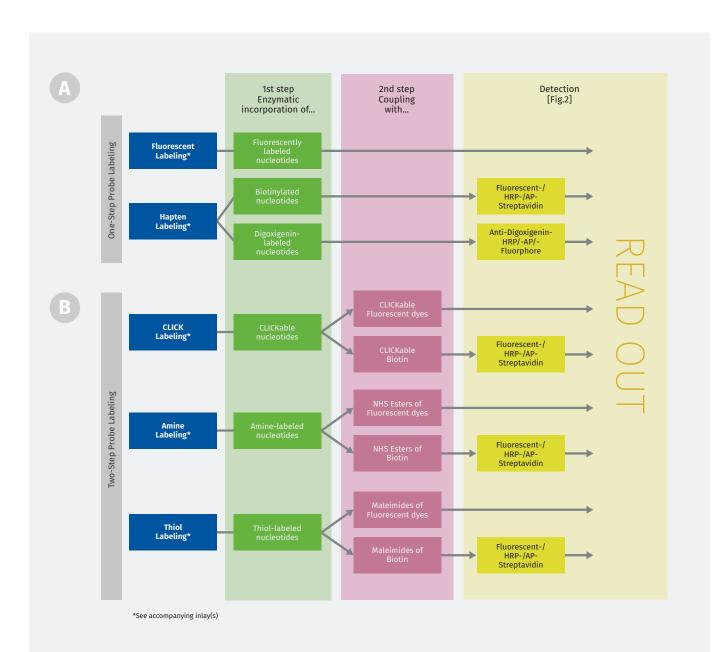
Figure 2

Fluorescently labeled probes can be detected directly after incorporation (A) whereas indirect detection via Biotin/Streptavidin (B.1) or Digoxigenin/Antibody (B.2) systems offers signal amplification and increased stability. Commonly used reporter enzymes include horse-radish peroxidase (HRP) and alkaline phosphatase (AP) that generate signals through an enzymatic reaction with chemiluminescent or chromogenic substrates.

## Incorporation of labels into DNA/RNA

The incorporation of fluorophores and haptens into DNA and RNA hybridization probes is achieved by enzymatic labeling techniques using modified nucleotides as substitutes for their natural counterparts (Fig. 3). In a one-step procedure, accordingly labeled nucleotides are incorporated for immediate detection (Fig. 3A). Alternatively, two-step labeling is achieved via i) incorporation of a reactive group followed by ii) coupling with the desired label (Fig. 3B).

To ensure optimal substrate properties of such modified nucleotides, their core structure is derived from natural analogs with the desired moiety attached via a linker to one of the available modifiable nucleotide positions (base, ribose, phosphate). Both, the position of linker attachment and the type of linker, are critical factors affecting substrate properties and resulting labeling efficiency [6].



**Figure 3** One-step probe labeling (A) is performed by incorporation of nucleotide analogs carrying a label (dye or hapten) while two-step probe labeling (B) requires in a first step introduction of a reactive group (amine, thiol, or "Clickable" moiety) followed by coupling with an appropriate label. NHS: N-hydroxy-succinimide, DBCO: Dibenzylcyclooctyne, HRP: horseradish peroxidase, AP: alkaline phosphatase, Cu: copper. Please note that this is a non-comprehensive illustration of existing labeling methods.

The labeling strategy depends on the size and type of nucleic acid to be labeled (DNA or RNA), the required sensitivity for the downstream application and the availability of appropriately labeled nucleotide analogs for enzymatic incorporation. For example, labeling with near-infrared dyes is, on the one hand, of great importance when working with biological samples since near IR-fluorescence is not disturbed by biological autofluorescence (haemoglobins or cytochromes). Nucleotides labeled with near IR-dyes however, on the other hand, often show insufficient labeling efficiencies due to sterical hindrance caused by

the bulky fluorophore. This problem can be circumvented by the use of smaller labels such as haptens or reactive group-carrying nucleotides followed by coupling with a near-IR dye (Fig. 3). Incorporation of modified nucleotides into DNA & RNA can be done random (i.e. incorporation of multiple nucleotides in a random fashion along the entire nucleic acid) or by selective end-labeling (incorporation of one or multiple nucleotides exclusively at the 3´- and/or 5´-end). Random labeling usually achieves higher labeling densities and thus higher sensitivities (Tab. 1).

#### Table 1 Standard enzymatic procedures are available for the preparation non-radioactively labeled probes.

Taq: Thermus aquaticus; Vent exo-: Thermococcus litoralis exo- DNA polymerase; MMLV: Moloney Murine Leukemia Virus; AMV: Avian Murine Virus; exo-: lack of exonuclease activity.

Template	Method	Labeled Probe	Labeling site	Enzyme	References
DNA	PCR	DNA	random	Thermophilic polymerases e.g. Taq Polymerase or Vent exo	[7], [8]
	Nick Translation	DNA	random	DNAse I/ DNA Polymerase I	[8], [9], [10]
	Primer Extension	DNA	random	Klenow Fragment 3'->5' exo <sup>-</sup> Taq Polymerase	[8]
	3´-End Labeling	DNA	3´-OH	Terminal Deoxynucleotidyl Transferase (TdT)	[11], [12]
	5´-End Labeling	DNA	5´-OH	T4 Polynucleotide Kinase (T4 PNK)	[13]
	In vitro Transcription	cRNA	random	SP6 RNA Polymerase T7 RNA Polymerase T3 RNA Polymerase	[14], [15]
	·	cRNA	5´-OH	T7 RNA Polymerase	[16]
RNA	Reverse Transcription	cDNA	random	MMLV Reverse Transcriptase AMV Reverse Transcriptase	[7], [14]
	3´-End Labeling	RNA	3´-OH	Terminal Deoxynucleotidyl Transferase (TdT)	[17], [18]
		RNA	3´-OH	Yeast Poly A Polymerase	[19]
		RNA	3´-OH	T4 RNA Ligase	[20]
	5´-End Labeling	RNA	5´-OH	T4 Polynucleotide Kinase (T4 PNK)	[13], [21]

#### Recommended reading:

- [1] Besse et al. (2009) Drosophila PTB promotes formation of high-order RNP particles and represses oskar translation. Genes Dev. 23:195.
- [2] Diamandis et al. (1991) The biotin-(strept)avidin system: principles and applications in biotechnology. Clin. Chem. 37:625.
- [3] McQuaid et al. (1995) A comparison of digoxigenin and biotin labelled DNA and RNA probes for in situ hybridization. Biotech Histochem. 70:147.
- [4] Chevalier et al. (1997) Biotin and digoxigenin as labels for light and electron microscopy in situ hybridization probes: where do we stand? J. Histochem. Cytochem. 45:481.
- [5] Leary et al. (1983) Rapid and sensitive colorimetric method for visualizing biotin-labeled DNA probes hybridized to DNA or RNA immobilized on nitrocellulose: Bio-blots. Proc. Natl. Acad. Sci. USA 80:4045
- [6] McGall (2005) Nucleoside triphosphate analogs for nonradioactive labeling of nucleic acids. In: Nucleoside Triphosphates and their Analogs (Vaghefi). CRC Press.
- [7] Anderson et al. (2005) Incorporation of reporter-labeled nucleotides by DNA polymerases. Biotechniques 38:257.
- [8] Yu et al. (1994) Cyanine dye dUTP analogs for enzymatic labeling of DNA probes. Nucleic Acids Res. 22:3226.
- [9] Langer et al. (1981) Enzymatic synthesis of biotin-labeled polynucleotides: novel nucleic acid affinity probes. Proc. Natl. Acad. Sci. USA 78:6633.
- [10] Gebeyehu et al. (1987) Novel biotinylated nucleotide-analogs for labeling and colorimetric detection of DNA. Nucleic Acids Res. 15:4513.
- [11] Schmitz et al. (1991) Nonradioactive labeling of oligonucleotides in vitro with the hapten digoxigenin by tailing with terminal transferase. Anal. Biochem. 192:222.
- [12] Flickinger et al. (1992) Differential incorporation of biotinylated nucleotides by terminal deoxynucleotidyl transferase. Nucleic Acids Res. 20:2382.
- [13] Grant et al. (2007) A facile method for attaching nitroxide spin labels at the 5' terminus of nucleic acids. Nucleic Acids Res. 35:e77.
- [14] Richter et al. (2002) Comparison of fluorescent tag DNA labeling methods used for expression analysis by DNA microarrays. Biotechniques 33:620.
- [15] t Hoen et al. (2003) Fluorescent labelling of cRNA for microarray applications. Nucleic Acids Res. 31:e20.
- [16] Samanta et al. (2014) A modified dinucleotide for site-specific RNA-labeling by transcription priming and click chemistry. Chem. Communication. 50:1313.
- [17] Zhao et al. (2012) Simple and sensitive microRNA labeling by terminal deoxynucleotidyl transferase. Acta Biochim. Biophys. Sin. (Shanqhai) 44:129.
- [18] Rosemeyer et al. (1995) Nonradioactive 3'-end-labeling of RNA molecules of different lengths by terminal deoxynucleotidyltransferase. Anal. Biochem. 224:446.
- [19] Martin et al. (1998) Tailing and 3'-end labeling of RNA with yeast poly(A) polymerase and various nucleotides. RNA 4:226.
- [20] Wang et al. (2007) Direct and sensitive miRNA profiling from low-input total RNA. RNA 13:151.
- [21] Akabayov et al. (2002) Immobilization of RNA: Application to Single Molecule Spectroscopy. Molecular, Cellular and Tissue Engineering. Proceedings of the IEEE-EMBS Special Topic Conference 71.



