

Jena Bioscience



Fluorescent Biomolecules and Their Building Blocks – Design and Applications

Fluorescent Nucleotides: A powerful toolbox for labeling of biological macromolecules

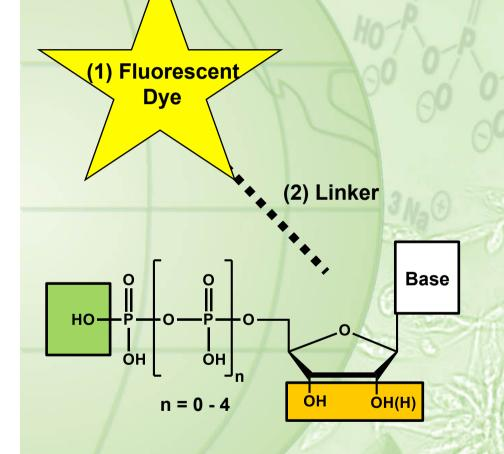
July 5th - 8th 2012

Jena Bioscience GmbH Loebstedter Str. 80 07749 Jena, Germany Tel.: +49-3641-628-5000 Fax: +49-3641-628-5100 e-Mail: info@jenabioscience.com

Many biological macromolecules are accessible via nucleotides

	DNA & RNA	 made of nucleotides (deoxy- or ribonucleotides) hybridize with complementary oligo probes
	Proteins /	Several thousands <i>interact</i> with nucleotides or DNA/ RNA:
	Enzymes	 Major signal transduction pathways require NTPs as cofactors GPCR/G-protein ←> GTP ←> 60% of current pharmaceutical drugs⁽¹⁾ Protein Kinases ←> ATP
		 Polymerases ← → dNTPs Motor proteins, chaperones, transcription factors, small molecule kinases,
	Fluorescent	nucleotides facilitate monitoring physiological processes

Conservative estimate: There are millions of fluorescent nucleotides possible



- (1) >100 fluorescent dyes commercially available
- (2) approx.10 linkers attachable to ~5 positions of nucleotide
- (3) 5 natural bases, each with >10 known modifications
- (4) 2 natural sugars, each with >10 known modifications
- (5) 3 natural phosphate states (mono-, di-, triphosphate) ignoring higher phosphates and substitutions of oxygen with sulfur or –NH– / – CH₂–

15 million structures of fluorescent nucleotides

Do we really need so many...?

No, we don't...we only need two general types of fluorescent nucleotide probes

A set of pre-designed probes available from shelf for common applications - presumably a few hundred...thousand are sufficient -

&

Tailor-made probes

for particular, special purposes



Site (position) of label attachment



Linkers: More than just a handle



Labeling reaction: Simply CLICK it!

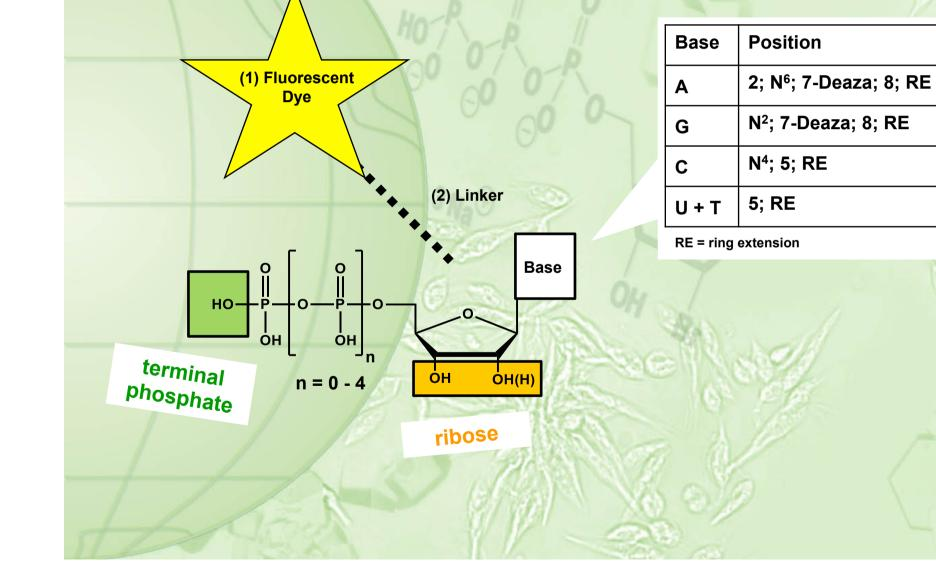


Selected applications



Label position: Base, sugar or terminal phosphate For base modification purines offer more options than pyrimidines

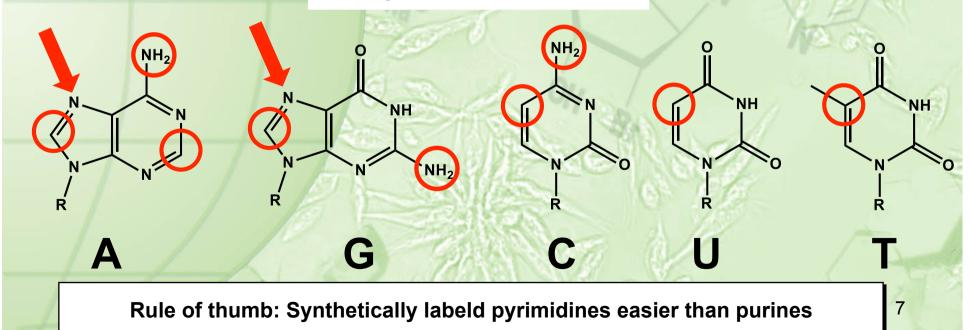
6



A has 4 labeling sites + ring extensions, G has 3 +1

Base	Position
Α	2; N ⁶ ; 7-Deaza; 8; RE
G	N²; 7-Deaza; 8; RE
С	N⁴; 5; RE
U + T	5; RE

RE = ring extension





Site (position) of label attachment



Linkers: More than just a handle

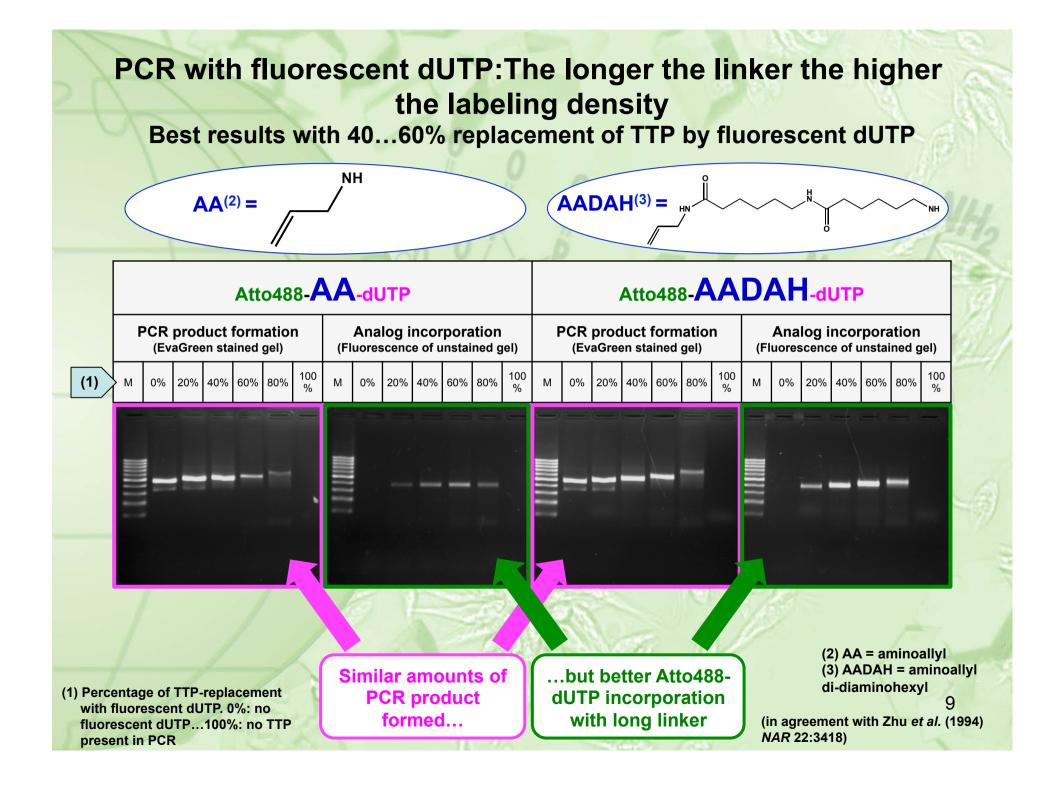


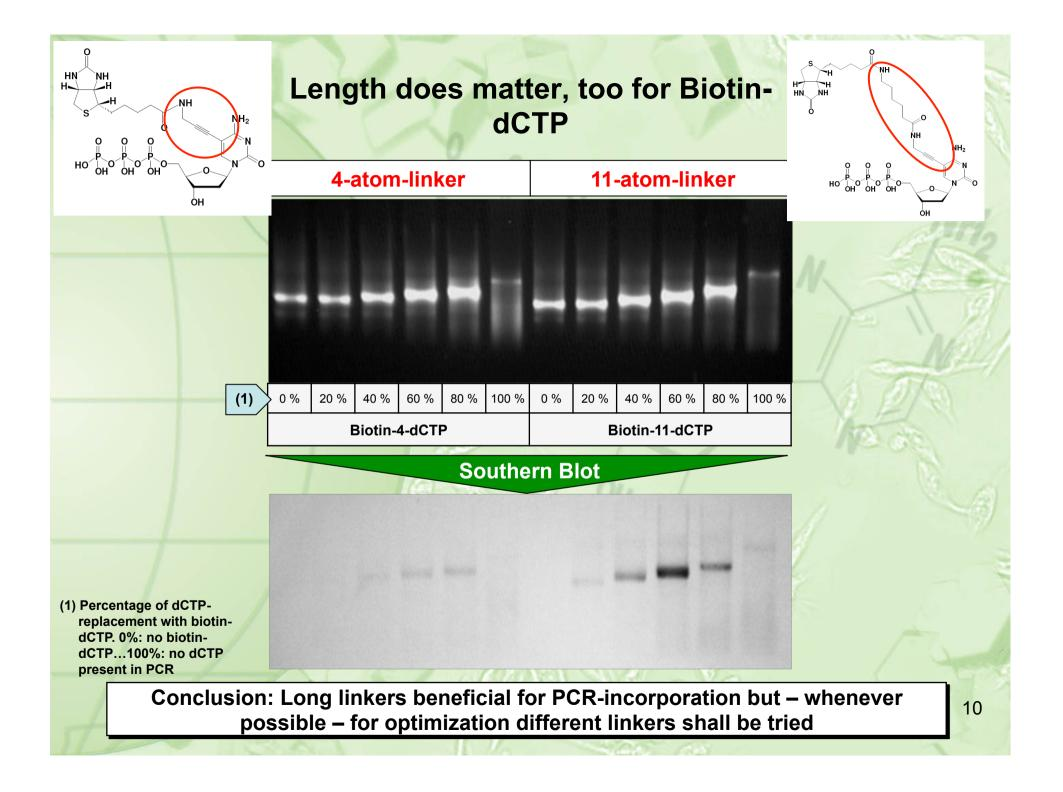
Labeling reaction: Simply CLICK it!



Selected applications







11



Site (position) of label attachment



Linkers: More than just a handle



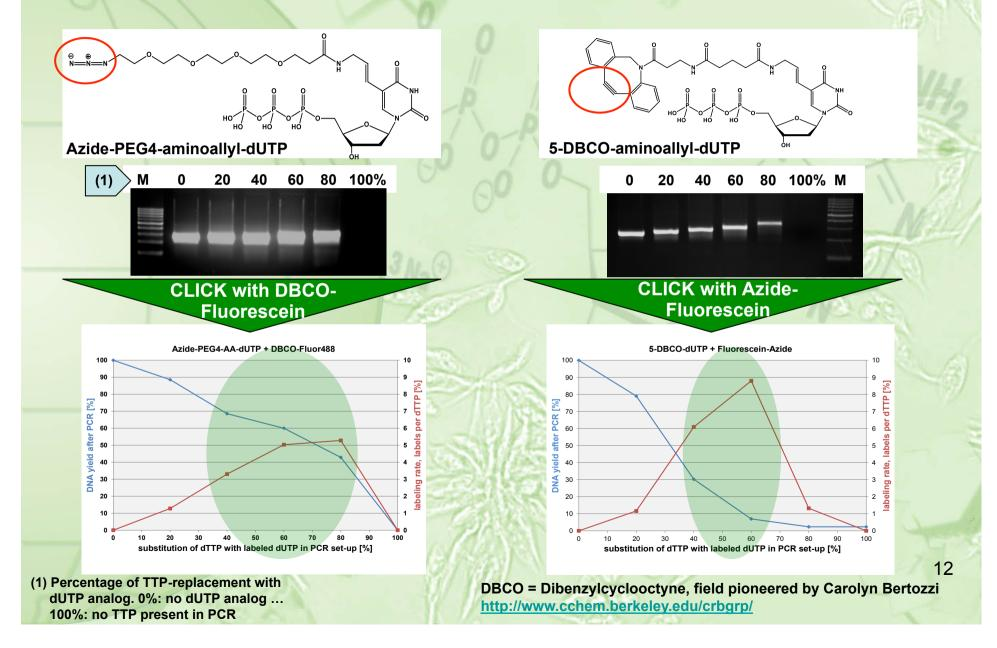
Labeling reaction: Simply CLICK it!



Selected applications



Classic Amine-NHS-ester chemistry can be replaced by CLICK reactions





Site (position) of label attachment



Linkers: More than just a handle



Labeling reaction: Simply CLICK it!

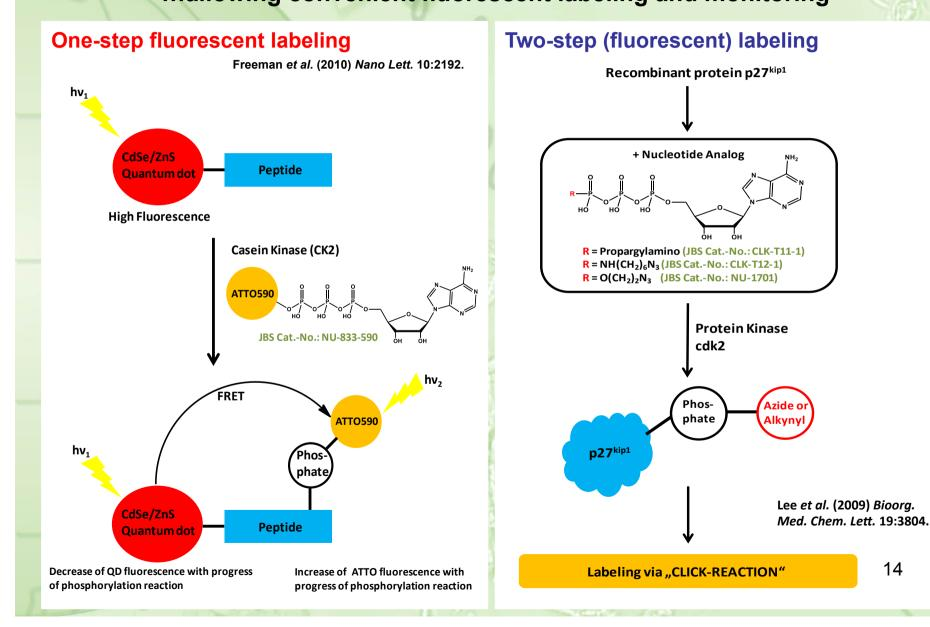


Selected applications



Protein kinases transfer modified y-phosphate of ATP to their substrates... ...allowing convenient fluorescent labeling and monitoring

14





Site (position) of label attachment



Linkers: More than just a handle



Labeling reaction: Simply CLICK it!



Selected applications



Nucleotides and their analogs is largest of eight lines of business



Roughly 2.000 nucleotides in stock & custom projects (lead discovery/optimization for big pharma but also academia-compatible custom syntheses) ¹⁶

Easy navigation by structure, application & structural formula



Structure Search beta

Search all our small molecule products (including all nucleosides and nucleotides) by sub-structure.

Nucleotides by Structure

Fluorescent Nucleotides

- Dye-/Quencher-labeled Nucleotides
- Intrinsically Fluorescent Analogs
- Nucleotide Probes Trove

Non-hydrolyzable Nucleotides

- <u>α-Phosphate modified Nucleotides</u>
- <u>α,β non-hydrolyzable Nucleotides</u>
- β-Phosphate modified Nucleotides
- β, y non-hydrolyzable Nucleotides
- y-Phosphate modified Nucleotides
- Non-hydrolyzable Di-Nucleotides
- Non-hydrolyzable Nucleotide Kits

Nucleotides labeled with...

- Biotin
- Digoxigenin
- Photo-labile groups ("Caged")
- Free amino group (-NH₂)
- Redox Dyes
- Halogen atoms (F, Cl, Br, I)
- Mercury (Hg)
- Selen (Se)

Analogs and Derivatives of...

- Cap
- Puromycin
- Coenzyme A (CoA)
- <u>NAD</u>

Unmodified (natural) Nucleotides

- <u>dNTPs</u>
- <u>dNDPs</u>
- NTPs
- NDPs
- <u>NMPs</u>

Important Structure Motifs

- Cyclic Nucleotides
- <u>Di-nucleoside Phosphates</u>
- 6-Thio Guanosines and Inosines
- <u>7-Deaza Purines</u>
- <u>7-Methyl Guanosines</u>
- <u>4-Substituted Pyrimidines</u>
- <u>5-Methyl Cytidines and related</u>
- Unmodified and Modified ddNTPs

Nucleotide Trove

- 6-Modified Purines
- <u>8-Oxo Guanosines</u>
- <u>2'-Deoxy Uridines</u>
- <u>3'-Deoxy Nucleotides</u>
- Nucleoside Bisphosphates
- ara-Nucleotides
- Unmodified Purines
- Modified dNTPs
- Miscellaneous

📆 Custom Synthesis

Didn't find the nucleotide analog you're looking for? Contact us at <u>nucleotides@jenabioscience.com</u> for a custom synthesis! With our pre-made building blocks and in-house expertise we manufacture even the most exotic nucleotide analog from the mg to kg scale - find more information in the PDF file above.

Nucleotides by Application

...on DNA

- PCR & Co.
- Sequencing & Genotyping
- Mutagenesis
- Crosslinking

...on RNA

- in vitro Synthesis
- <u>5'-Capping</u>

...on Proteins & Enzymes

- <u>Co-translational Protein Labeling</u>
- Kinase Signaling
- G-Protein Signaling
- Tubulin Assembling
- Structural Biology

... in Cell Biology

- <u>Cell Cycle & Proliferation</u>
- Apoptosis
- Protein-DNA/-RNA Interaction (EMSA)
- Epigenetics

... in Drug Discovery

- <u>SELEX/Aptamer Modification</u>
- mRNA Modification
- Antiviral Nucleotides
- Cytostatic Nucleotides





Navigation



Structure Search^{beta} Search all our small molecule products (including all nucleosides and nucleotides) by

Nucleotides by Structure

Fluorescent Nucleotides

- Dye-/Quencher-labeled Nucleotides
- Intrinsically Fluorescent Analogs
- Nucleotide Probes Trove

Non-hydrolyzable Nucleotides

- α-Phosphate modified Nucleotides
- α,β non-hydrolyzable Nucleotides
- β-Phosphate modified Nucleotides
- β, y non-hydrolyzable Nucleotides
- y-Phosphate modified Nucleotides
- Non-hydrolyzable Di-Nucleotides
- Non-hydrolyzable Nucleotide Kits

Nucleotides labeled with...

- Biotin
- Digoxigenin
- Photo-labile groups ("Caged")
- Free amino group (-NH₂)
- Redox Dyes
- Halogen atoms (F, Cl, Br, I)
- Mercury (Hg)
- Selen (Se)

Analogs and Derivatives of ...

- Cap
- Puromycin
- Coenzyme A (CoA)
- NAD

Unmodified (natural) Nucleotides

- dNTPs
- dNDPs
- NTPs
- NDPs

📆 Custom Synthesis

Didn't find the nucleotide analog you're looking for? Contact us at <u>nucleotides@jenabioscience.com</u> for a custom synthesis! With our pre-made building blocks and in-house expertise we manufacture even the most exotic nucleotide analog from the mg to kg scale - find more information in the PDF file above.

Nucleotides by Application

...on DNA

- PCR & Co.
- Sequencing & Genotyping
- Mutagenesis
- Crosslinking

...on RNA

- in vitro Synthesis
- <u>5'-Capping</u>

...on Proteins & Enzymes

- <u>Co-translational Protein Labeling</u>
- Kinase Signaling
- G-Protein Signaling
- Tubulin Assembling
- Structural Biology

... in Cell Biology

- Cell Cycle & Proliferation
- Apoptosis
- Protein-DNA/-RNA Interaction (EMSA)
- Epigenetics

... in Drug Discovery

- SELEX/Aptamer Modification
- mRNA Modification
- Antiviral Nucleotides
- Cytostatic Nucleotides

20