



Fluorescent and Hapten-Labeled Nucleotide Analogs

Labeled nucleotides are critical elements for sequence detection in a wide variety of techniques including *in situ* hybridization, microarrays and DNA sequencing. Our proprietary synthesis and purification delivers consistently high purity products to ensure the most efficient incorporation for optimum results in your experiment. We offer a broad array of standard products as well as custom synthesis and packaging options.

Labeled Nucleotides for Direct Fluorescence Detection

Dideoxynucleotides

>95% pure by analytical HPLC.

Fluorescein-Labeled: 1 mM, 25 nmol

Cat. No.	Product Description
NEL400001EA	Fluorescein-12-ddCTP
NEL401001EA	Fluorescein-12-ddUTP
NEL402001EA	Fluorescein-12-ddATP

ROX-Labeled: 0.1 mM, 2.5 nmol

Cat. No.	Product Description
NEL476001EA	ROX-ddUTP
NEL478001EA	ROX-ddATP

R6G-Labeled: 0.1 mM, 2.5 nmol

Cat. No.	Product Description
NEL489001EA	R6G-ddCTP
NEL490001EA	R6G-ddATP

TAMRA-Labeled: 0.1 mM, 2.5 nmol

Cat. No.	Product Description
NEL472001EA	TAMRA-ddUTP
NEL473001EA	TAMRA-ddCTP
NEL474001EA	TAMRA-ddATP
NEL475001EA	TAMRA-ddGTP

Cyanine 3-Labeled: 0.1 mM, 2.5 nmol

Cat. No.	Product Description
NEL586001EA	Cyanine 3-ddATP
NEL587001EA	Cyanine 3-ddGTP

Cyanine 5-Labeled: 0.1 mM, 2.5 nmol

Cat. No.	Product Description
NEL588001EA	Cyanine 5-ddCTP
NEL589001EA	Cyanine 5-ddUTP
NEL591001EA	Cyanine 5-ddGTP

R110-Labeled: 0.1 mM, 2.5 nmol

Cat. No.	Product Description
NEL492001EA	R110-ddUTP
NEL493001EA	R110-ddCTP
NEL495001EA	R110-ddGTP

Deoxynucleotides

>95% pure by analytical HPLC. 1 mM, 25 nmol package sizes, unless otherwise noted.

dUTP Analogs

Cat. No.	Product Description
NEL413001EA	Fluorescein-12-dUTP
NEL417001EA	Texas Red®-5-dUTP
NEL418001EA	Lissamine-5-dUTP
NEL455001EA	Diethylaminocoumarin-5-dUTP
NEL578001EA	Cyanine 3-dUTP
NEL579001EA	Cyanine 5-dUTP
NEL621001EA	Smart Pack dUTP (5 x 25 nmol Cyanine 3-dUTP and 5 x 25 nmol Cyanine 5-dUTP)

dCTP Analogs

Cat. No.	Product Description
NEL424001EA	Fluorescein-12-dCTP
NEL426001EA	Texas Red®-5-dCTP
NEL576001EA	Cyanine 3-dCTP
NEL577001EA	Cyanine 5-dCTP
NEL620001EA	Smart Pack dCTP (5 x 25 nmol Cyanine 3-dCTP and 5 x 25 nmol Cyanine 5-dCTP)

dATP Analogs

Cat. No.	Product Description
NEL465001EA	Fluorescein-12-dATP
NEL471001EA	Texas Red®-5-dATP
NEL592001EA	Cyanine 3-dATP
NEL593001EA	Cyanine 5-dATP

dGTP Analogs

Cat. No.	Product Description
NEL429001EA	Fluorescein-12-dGTP
NEL594001EA	Cyanine 3-dGTP
NEL595001EA	Cyanine 5-dGTP

Ribonucleotides

>95% pure by analytical HPLC. 1 mM, 25 nmol package sizes, unless otherwise noted.

UTP Analogs

Cat. No.	Product Description
NEL414001EA	Fluorescein-12-UTP*
NEL582001EA	Cyanine 3-UTP*
NEL583001EA	Cyanine 5-UTP*
NEL623001EA	Smart Pack UTP* (5 x 100 nmol Cyanine 3-UTP and 5 x 100 nmol Cyanine 5-UTP)

*Available as: 10 mM, 100 nmol package

CTP Analogs

Cat. No.	Product Description
NEL434001EA	Fluorescein-12-CTP*
NEL580001EA	Cyanine 3-CTP*
NEL581001EA	Cyanine 5-CTP*

*Available as: 10 mM, 100 nmol package

ATP Analogs

Cat. No.	Product Description
NEL439001EA	Fluorescein-12-ATP

GTP Analogs

Cat. No.	Product Description
NEL496001EA	Fluorescein-12-GTP

Labeled Nucleotides for Indirect Detection Methods

Deoxynucleotides hapten-labeled: 1 mM, 25 nmol, >95% pure by analytical HPLC

Cat. No.	Product Description
NEL538001EA	Biotin-11-dCTP
NEL539001EA	Biotin-11-dUTP
NEL540001EA	Biotin-11-dATP
NEL541001EA	Biotin-11-dGTP
NEL551001EA	DNP-11-dUTP
NEL562001EA	Aminodigoxigenin-9-dCTP

Dideoxynucleotides hapten-labeled: 1 mM, 25 nmol, >95% pure by analytical HPLC

Cat. No.	Product Description
NEL546001EA	Biotin-11-ddCTP
NEL547001EA	Biotin-11-ddUTP
NEL548001EA	Biotin-11-ddATP
NEL549001EA	Biotin-11-ddGTP
NEL573001EA	Aminodigoxigenin-9-ddGTP

Ribonucleotides hapten-labeled: 1 mM, 25 nmol >95% pure by analytical HPLC

Cat. No.	Product Description
NEL542001EA	Biotin-11-CTP
NEL543001EA	Biotin-11-UTP
NEL544001EA	Biotin-11-ATP
NEL545001EA	Biotin-11-GTP
NEL555001EA	DNP-11-UTP

Complementary Products

Cat. No.	Product Description
NEF709001EA	Antifluorescein-AP
NEF710001EA	Antifluorescein-HRP
NEL103001EA	Western Lightning Plus Chemiluminescent HRP Substrate
NEL300001EA	4CN Plus Chromogenic HRP Substrate
NEL601001KT	CDP-Star [®] Chemiluminescent AP Substrate
NEL720001EA	Streptavidin-Fluorescein
NEL721001EA	Streptavidin-Texas Red [®]
NEL722001EA	Streptavidin-Coumarin
NEL750001EA	Streptavidin-HRP
NEL751001EA	Streptavidin-AP
NEL937001EA	BCIP/NBT-AP Substrate
NEL938001EA	DAB-HRP Substrate

Excitation and Emission Spectra for Common Fluorophores

Fluorophore	Excitation (nm)	Emission (nm)
Pyrene	340	376, 395
Coumarin	402	443
Diethylaminocoumarin	426	480
FAM	490	520
Fluorescein Chlorotriazinyl	492	514
Fluorescein	494	517
R110	500	525
Eosin	517	537
JOE	524	550

Fluorophore	Excitation (nm)	Emission (nm)
R6G	525	550
Cyanine 3	550	568
Tetramethylrhodamine	550	570
TAMRA	552	575
Lissamine [®]	570	588
ROX	580	605
Texas Red [®]	593	612
Naphthofluorescein	590	667
Cyanine 5	650	668

OnPoint Custom Services

PerkinElmer will custom synthesize and purify labeled nucleotides for your research. Our proprietary purification system provides analogs of the highest purity available. In many cases, the analog can be provided isomerically pure.

PerkinElmer can accommodate your nucleotide needs. Larger package sizes may be arranged for any product listed. Buffers and analog concentrations other than described are also available.

For more information on PerkinElmer Fluorescent and Hapten-labeled nucleotide analogs, visit

www.perkinelmer.com/category/fluorescent-hapten-nucleotides

This product or the use of this product may be covered by one or more patents owned by PerkinElmer LAS, Inc. including U.S. Patent Nos. 5,242,796; 5,306,618; 5,332,666; 5,558,991; 5,608,063; and 5,625,081, or foreign equivalents thereof. Those products incorporating a cyanine dye may be additionally covered by one or more of the following issued US patents owned by PerkinElmer LAS, Inc. including US Patent Nos. 6,114,350; 6,197,956; 6,204,389; 6,224,644; and 6,437,141 or foreign equivalents thereof on the cyanine dye precursors, uses and labeled moieties. PerkinElmer makes no warranty that this product may be used in any way without infringing the patent or other intellectual property rights of a third party.

PerkinElmer, Inc.
940 Winter Street
Waltham, MA 02451 USA
P: (800) 762-4000 or
(+1) 203-925-4602
www.perkinelmer.com



For a complete listing of our global offices, visit www.perkinelmer.com/ContactUs

Copyright © 2008-2018, PerkinElmer, Inc. All rights reserved. PerkinElmer[®] is a registered trademark of PerkinElmer, Inc. All other trademarks are the property of their respective owners.

007384D_01 PKI