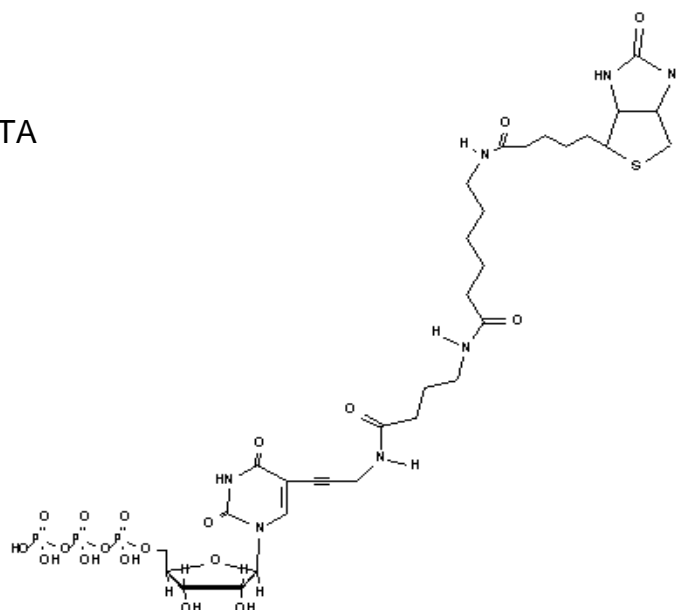


Caution: For Laboratory Use. A product for research purposes only.

BIOTIN-16-UTP

Product Number: NEL 630

QUANTITY: 250 nmol
FORM: 25 μ L solution
CONCENTRATION: 10.0 mM
SOLVENT: 10 mM Tris-HCl, pH 7.6, 1 mM EDTA
EXTINCTION COEFFICIENT: 13,000 M⁻¹cm⁻¹
 (289 nm, Phosphate buffer, pH 7)



EXCITATION MAXIMUM: 289 nm

INTRODUCTION

Nucleotide analogs^{1,2,3} are biologically active with a variety of DNA and/or RNA polymerases. Labeling methods such as: nick translation, random priming, polymerase chain reaction, 3'-end labeling, or transcription of RNA using SP6, T3, or T7 RNA polymerases may be used. Some analogs demonstrate variations in relative performance depending upon nucleotide and label (fluorophore or hapten) selected due to enzyme preferences. Labeled probes may be used in applications including (but not limited to) chromosome mapping. These analogs are intended to be detected either directly by their fluorescence when using a fluorescently labeled analog or indirectly when appropriately labeled antibodies or streptavidin are available. Indirect detection may be either colorimetric, chemiluminescence, or fluorescence. Signal amplification may be obtained using NEN's patented Tyramide

Signal Amplification process (TSA™). **For additional information: call 1-800-762-4000 or visit our WEB site at <http://las.perkinelmer.com>.**

QUALITY CONTROL

The analog is purified by HPLC chromatography. Analytical HPLC is done to ensure initial purity is >95%. UV/VIS absorption spectra are obtained in aqueous phosphate buffer and used to determine concentration. Copies of representative spectra, labeling protocols, and information about TSA™ are available from Technical Service at 1-800-551-2121 or visit our web site: <http://www.perkinelmer.com>.

STABILITY AND STORAGE CONDITIONS

Nucleotides labeled with fluorophores should be protected from extended exposure to light. These nucleotide analogs are stable kept in a refrigerator or colder for at least 1 year. Minimizing freeze-thaw cycles and exposure to light are the most critical factors to consider for long term usage.

For Research Use Only:

¹Those products incorporating a cyanine dye are covered under the following issued US Patent Nos: 6114350, 6197956, 6204389, and 6224644 on the cyanine dye precursors, uses and labeled moieties.

²This product may not be used for DNA sequencing unless (a) used with a DNA sequencer instrument purchased from PerkinElmer Health Sciences, Inc. or its sublicensees, or (b) a separate license for such use is obtained from Applied Biosystems, Inc., Foster City, CA.

³The use of this product for primer extension may be covered by one or more of the following US patents (or their foreign counterparts): 5888819, 5952174, 6004744, and/or 6013431, and to the extent covered may not be used unless a separate license for such use is obtained from Beckman Coulter, Inc., of Fullerton, CA.

PerkinElmer, Inc.
549 Albany Street
Boston, MA 02118 USA
P: (800) 762-4000 or (+1) 203-925-4602
www.perkinelmer.com/nenradiochemicals

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

Copyright ©2010, PerkinElmer, Inc. All rights reserved. PerkinElmer® is a registered trademark of PerkinElmer, Inc. All other trademarks are the property of their respective owners.

