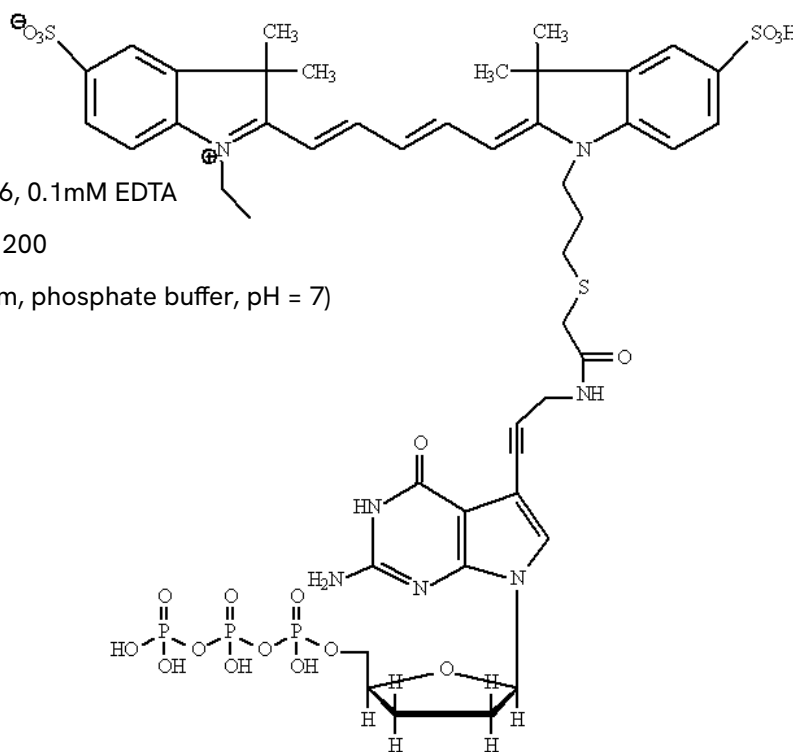


Cyanine 5-ddGTP

Product Number: NEL591001EA

Quantity: 2.5 nmol
 Form: 25 μ L solution
 Concentration: 0.1 mM
 Solvent: 10 mM Tris-HCL, pH 7.6, 0.1mM EDTA
 Formula: $C_{46}H_{56}N_7O_{19}P_3S_3$. MW 1200
 Extinction Coefficient: 250,000 $M^{-1}cm^{-1}$ (650nm, phosphate buffer, pH = 7)
 Excitation Maximum: 650 nm
 Emission Maximum: 668 nm



Introduction

Fluorescent dideoxynucleotide analogs are 3'-end chain terminators used for DNA sequencing, RFLP mapping, DNA fingerprinting, and point mutation screening analysis. Labeled DNA patterns may be obtained by either separating labeled fragments using standard polyacrylamide gel electrophoresis techniques or with primer/template coated microarrays using 96-well microplates or glass slides as solid supports. Detection is via the direct fluorescence of the nucleotide analog using classical excitation/emission or when using two fluorescent dyes in combination via fluorescence resonance energy transfer (FRET). These analogs are available with a variety of fluorophores attached to each of the four dideoxynucleotides to permit maximum assay flexibility.

Quality Control

The nucleotide analog is purified by HPLC chromatography. Analytical HPLC is used as a quality control check to ensure chemical and isomeric purity >95%. UV/VIS absorption spectra are obtained in aqueous phosphate buffer to determine concentration. Relative fluorescence quantum yields are not necessarily the same for the four different base nucleotide analogs.

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 most critical factors to consider for long term usage.

The information provided in this document is for reference purposes only and may not be all-inclusive. Revvity, Inc., its subsidiaries, and/or affiliates (collectively, "Revvity") do not assume liability for the accuracy or completeness of the information contained herein. Users should exercise caution when handling materials as they may present unknown hazards. Revvity shall not be liable for any damages or losses resulting from handling or contact with the product, as Revvity cannot control actual methods, volumes, or conditions of use. Users are responsible for ensuring the product's suitability for their specific application. REVVITY EXPRESSLY DISCLAIMS ALL WARRANTIES, INCLUDING WARRANTIES OF MERCHANTABILITY OR FITNESS FOR A PARTICULAR PURPOSE, REGARDLESS OF WHETHER ORAL OR WRITTEN, EXPRESS OR IMPLIED, ALLEGEDLY ARISING FROM ANY USAGE OF ANY TRADE OR ANY COURSE OF DEALING, IN CONNECTION WITH THE USE OF INFORMATION CONTAINED HEREIN OR THE PRODUCT ITSELF

NEL591001EA-R REV01

The logo for Revvity, featuring the word "revvity" in a lowercase, sans-serif font.