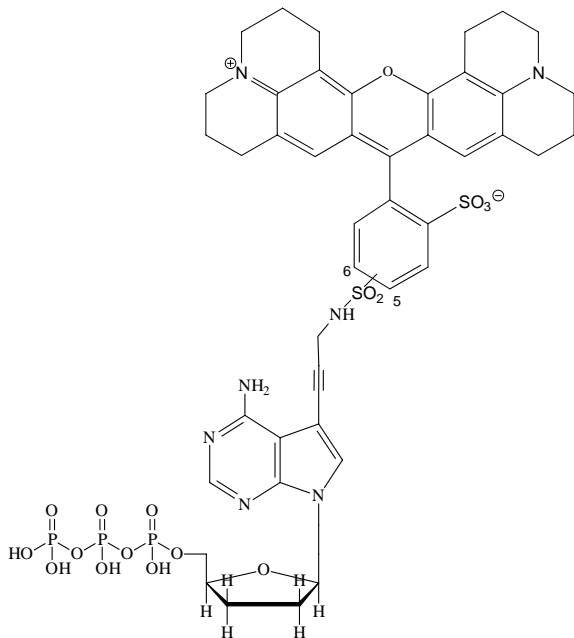
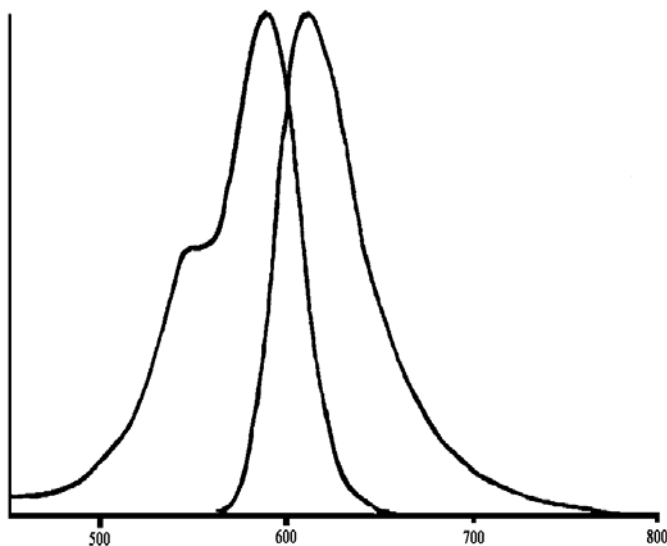


**Texas Red®-5-ddATP
NEL411**



QUANTITY: 25 nmol
FORM: 25 µL solution
CONCENTRATION: 1.0 mM
SOLVENT: 10 mM Tris-HCl, pH 7.6, 1 mM EDTA
FORMULA: C₄₆H₅₁N₇O₁₇P₃S₂ FW = 1131
COLOR: bluish red
EXTINCTION COEFFICIENT: 85,000 M⁻¹cm⁻¹
 (593 nm, Phosphate buffer, pH = 7)



WAVELENGTH: EXCITATION 593nm
 (Maxima) EMISSION 612nm

PC2418 – 0105

INTRODUCTION

Fluorescent 2',3'-dideoxynucleotide analogs^{1,2} are biologically active with a variety of DNA polymerases. Since dideoxynucleotides are chain terminators, they result in the analog being extended in a base specific manner onto the 3'-end of the DNA chain. Dideoxynucleotide analogs may be used in a variety of applications which allow determination of a genetic profile based on single nucleotide polymorphisms (SNP)². Analogs labeled with a fluorophore are intended to be detected directly by their fluorescence properties. **For additional information: call 1-800-762-4000 or visit our WEB site at http://www.perkinelmer.com/nucleotide_analogs.**

QUALITY CONTROL

The nucleotide analog is purified by HPLC chromatography. Analytical HPLC is used as a quality control check to ensure chemical 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.

For Research Use Only:

¹This 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. 5,047,519; 5,151,507; 5,558,991, and 5,608,063. Those products incorporating a cyanine dye are covered under the following issued US Patent Nos. 114,350, 6197, 956, 6,204,389, and 6,224,644 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 LAS, 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) – 5,888,819, 5,952,174, 6,004,744, 6,013,431 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.