



## 5-FAM-X-dUTP, lyophilized powder

Catalog Number	Packaging Size
C406B	25 nmol

**Storage upon receipt:** -20°C, protected from light.

### User Manual

GeneCopoeia, Inc.

9620 Medical Center Drive, #101

Rockville, MD 20850

USA

Tel: 1-866-360-9531; 1-301-762-0888

Fax: 1-301-762-3888

Inquiry: [inquiry@genecopoeia.com](mailto:inquiry@genecopoeia.com)

Order: [orders@genecopoeia.com](mailto:orders@genecopoeia.com)

Web: [www.genecopoeia.com](http://www.genecopoeia.com)

## I. Product Description

5-FAM-X-dUTP is supplied as lyophilized powder. The nucleotide is designed for enzymatic non-radioactive labeling of DNA by PCR, nick-translation, cDNA synthesis, random primed labeling, or primer extension. 5-FAM-X-dUTP can be enzymatically incorporated into DNA with Reverse Transcriptases, *Taq* DNA Polymerase, phi29 DNA Polymerase, Klenow Fragment, *exo-*, Klenow Fragment, and DNA Polymerase I. The X linker between FAM and dUTP can improve the incorporation efficiency.

## II. Specifications

Excitation/Emission (nm): 495/518 nm

Concentration: NA

Storage: Store at -20°C

Molecular Formula:  $C_{39}H_{38}N_4Na_3O_{21}P_3$

Molecular Weight: 1060.62

## III. General Characteristics

$\lambda_{max}=495\text{ nm}$ ,  $\epsilon=70.0 \times 10^3\text{ M}^{-1}\text{cm}^{-1}$  (pH 8.0).

