



## Andy Fluor 568-X-dUTP, 1 mM in TE buffer

Catalog Number	Packaging Size
C413A	25 µL

**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

Andy Fluor 568-X-dUTP is supplied as 1 mM solution in TE buffer. The nucleotide is designed for enzymatic non-radioactive labeling of DNA by PCR, nick-translation, cDNA synthesis, random primed labeling, or primer extension. Andy Fluor 568-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 Andy Fluor 568 and dUTP can improve the incorporation efficiency.

## II. Specifications

Excitation/Emission (nm): 578/602 nm

Concentration: 1 mM in TE buffer

Storage: Store at -20°C

## III. General Characteristics

$\lambda_{max}=578 \text{ nm}$ ,  $\epsilon=91.0 \times 10^3 \text{ M}^{-1}\text{cm}^{-1}$ .