

www.biotium.com

Revised: October 2, 2014

# **Product Information**

## Fluorescein-12-dUTP, lyophilized powder

Catalog Number 40063

Unit Size 25 nmole Molecular Weight ~991

Color and Form Orange solid

Absorption/Emission 494/521 nm

#### Storage and Handling

Store Fluorescein-12-dUTP at -20°C or -70°C, protected from light. When stored as recommended, product is stable for at least 6 months from date of receipt.

#### **Product Application**

Fluorescent conjugates of dUTP can be used for detection of apoptotic cells by direct fluorescence TUNEL labeling of DNA strand breaks in cells<sup>1</sup>, and can be used in place of dTTP in standard DNA labeling and synthesis protocols to generate fluorescent dsDNA and oligonucleotide probes.

### References

- 1. Gold et al. (1994). Lab Invest. 71 (2):219-25.
- 2. Slupphaug et al. (1993). Anal Biochem. 211 (1):164-9.
- 3. Hogrefe et al. (2002). PNAS 99 (2): 596-601.

Please visit our website at www.biotium.com to view our full selection of CF<sup>TM</sup> dye bioconjugates, including antibodies, antibody labeling kits, phalloidin, Annexin V and  $\alpha$ -bungarotoxin, as well as classic fluorescent nucleic acid dyes and antifade fluorescence mounting media.