

Product Information

Cyanine 555-dUTP, Lyophilized

Catalog Number 40064

Unit Size 25 nmole

Molecular Weight ~1131

Color and Form Red Solid

Absorption/Emission 550/570 nm

Storage and Handling

Store Cyanine 555-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 cells¹, 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™ dye bioconjugates, including antibodies, antibody labeling kits, phalloidin, annexin V and a-bungarotoxin, as well as classic fluorescent nucleic acid dyes and antifade fluorescence mounting media.