

Product Information

CF™555-ddCTP

Catalog Number: 40031

Unit Size: 25 nmol

Molecular Weight: ~1383

Color and Form: Red solid

Spectral Properties

$\lambda_{abs}/\lambda_{em} = 555/565$ nm (See Figure 1)

CF™555 is spectrally similar to Alexa Fluor®555, Cy®3, DyLight®549, and TRITC

Storage and Handling

Store desiccated at $\leq -20^{\circ}\text{C}$. When stored as directed, product should be stable for at least 6 months from the date of receipt. For aqueous solutions, prepare single use aliquots and store protected from light at -20°C for up to 6 months. Avoid freeze-thaw cycles. We recommend preparing a 1 mM stock solution in 10 mM Tris pH 7.4.

Product Application

Fluorophore conjugates of ddCTP can be used as 3'-end chain terminator in Sanger sequencing. CF555 has similar brightness and photostability as Alexa Fluor 555.

Related Products

Cat. #	Product description
40027	CF™555-dCTP, 25 nmol
40028	CF™647-dCTP, 25 nmol
40008	CF™488A-dUTP, 25 nmol
40002	CF™543-dUTP, 25 nmol
40005	CF™568-dUTP, 25 nmol
40006	CF™594-dUTP, 25 nmol
40007	CF™640R-dUTP, 25 nmol
40003	CF™680R-dUTP, 25 nmol

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 α -bungarotoxin, as well as reagents and kits for genomics and cell biology research.

CF™ dye technology is covered by pending US and international patents. Alexa Fluor is a registered trademark of Invitrogen; Cy is a registered trademark of GE Healthcare; DyLight is a registered trademark of Pierce Biotechnology.

Materials from Biotium are sold for research use only, and are not intended for food, drug, household, or cosmetic use.

Figure 1. Absorption/emission spectra of CF™555 conjugates

