

Product Information

CF™ Dye dCTP Conjugates

Catalog number	Size	Product	Molecular weight	Ex/Em (nm)
40067-T	5 nmol	CF™488A-dCTP	~1428	490/515
40067	25 nmol			
40057-T	5 nmol	CF™532-dCTP	~1447	527/558
40057	25 nmol			
40058-T	5 nmol	CF™543-dCTP	~1647	541/560
40058	25 nmol			
40027-T	5 nmol	CF™555-dCTP	~1553	555/565
40027	25 nmol			
40055-T	5 nmol	CF™568-dCTP	~1652	562/583
40055	25 nmol			
40056-T	5 nmol	CF™594-dCTP	~1667	593/614
40056	25 nmol			
40066-T	5 nmol	CF™640R-dCTP	~1593	642/662
40066	25 nmol			
40028-T	5 nmol	CF™647-dCTP	~1579	650/665
40028	25 nmol			
40068-T	5 nmol	CF™660R-dCTP	~1826	663/682
40068	25 nmol			

Storage and Handling

Store desiccated at ≤ -20°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

CF™ dyes are Biotium's next-generation fluorescent dyes, with combined advantages in brightness, photostability, and water solubility compared to other dyes like Alexa Fluor®, DyLight®, Cy® Dye, and IRDye®. CF dye conjugates of dCTP can be used in standard DNA labeling and synthesis protocols to generate fluorescent dsDNA and oligonucleotide probes.

Related Products

Catalog number	Product
30063	CF™488A TUNEL Assay Apoptosis Detection Kit
30064	CF™594 TUNEL Assay Apoptosis Detection Kit
30074	CF™640R TUNEL Assay Apoptosis Detection Kit
40004	CF™405S-dUTP
40008	CF™488A-dUTP
40002	CF™543-dUTP
40005	CF™568-dUTP
40006	CF™594-dUTP
40007	CF™640R-dUTP
40003	CF™680R-dUTP
40031	CF™555 ddCTP
40032	CF640R UTP
40029	Biotin-11-dUTP
40022	Biotin-16-dUTP
40030	Biotin-20-dUTP
40035	Biotin-11-CTP
40036	Biotin-11-dCTP
40033	Biotin-11-UTP
40023	Biotin-16-UTP
40034	Biotin-20-UTP
40020	5-Aminoallyl-dUTP
40021	5-Aminoallyl-UTP

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 fluorescent reagents and kits for genomics and cell biology research.

CF dye technology is covered by pending US and international patents. Alexa Fluor and DyLight are registered trademarks of Thermo Fisher Scientific; Cy Dye is a registered trademark of GE Healthcare; IRDye is a registered trademark of LI-COR Bioscience.

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