

Product Information

CF™ 640R-dCTP

Catalog Number: 40066

Unit Size: 25 nmol

Molecular Weight: ~1593

Color and Form: Blue solid

Spectral Properties

$\lambda_{abs}/\lambda_{em}$ = 642/662 nm (See Figure 1)

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 dCTP can be used in standard DNA labeling and synthesis protocols to generate fluorescent dsDNA and oligonucleotide probes.

Related Products

Cat. #	Product description
40027	CF™555-dCTP, 25 nmol
40028	CF™647-dCTP, 25 nmol
40031	CF™555-ddCTP, 25 nmol
40032	CF™640R-UTP, 25 nmol
40056	CF™594-dCTP, 25 nmol
40057	CF™532-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

Absorption/Emission Spectra of CF™640R Conjugated Antibodies

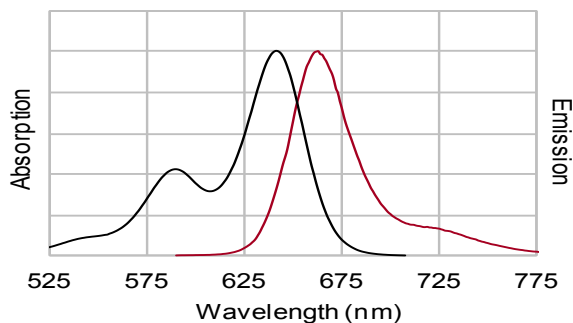


Figure 1. Absorption and emission spectra of CF™640R conjugated to goat anti-mouse IgG in PBS.

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.

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