

dCTP CF® Dye Conjugates

CF® dye dCTP can be used to synthesize labeled DNA probes for in-situ hybridization, microarray or blotting applications.



Product attributes

Colors Green, Orange, Red, Far-red

Product Description

CF® Dye dCTP can be used to synthesize labeled DNA probes for in situ hybridization, microarray, or blotting applications.

- Synthesize exceptionally bright and photostable fluorescent DNA probes
- Detect by fluorescence microscopy, flow cytometry, or in gel
- Compatible with fixed cells or tissue sections
- Choice of 9 CF® dye colors from green to far-red
- Biotium's CF® dyes have superior brightness, photostability, and are water-soluble

Supplied in lyophilized form. Also see our CF® Dye [dUTP Conjugates](#) and [TUNEL assay kits](#), available with a wide selection of bright and photostable CF® dyes.

Superior CF® Dyes

Biotium's next-generation CF® dyes were designed to be highly water-soluble with advantages in brightness and photostability compared to Alexa Fluor®, DyLight®, and other fluorescent dyes. Learn more about [CF® Dyes](#).

dCTP CF® Dye Conjugates

Product	Dye	Ex/Em	Size	SKU	Purchase
CF®488A dCTP 25 nmol	CF®488A 40067	490/515 nm Purchase 40067	5 nmol	40067-T	Purchase 40067-T
CF®532 dCTP 25 nmol	CF®532 40057	527/558 nm Purchase 40057	5 nmol	40057-T	Purchase 40057-T
CF®543 dCTP 25 nmol	CF®543 40058	541/560 nm Purchase 40058	5 nmol	40058-T	Purchase 40058-T
CF®555 dCTP 25 nmol	CF®555 40027	555/565 nm Purchase 40027	5 nmol	40027-T	Purchase 40027-T
CF®568 dCTP 25 nmol	CF®568 40055	562/583 nm Purchase 40055	5 nmol	40055-T	Purchase 40055-T
CF®594 dCTP ¹ 25 nmol	CF®594 40056	593/614 nm Purchase 40056	5 nmol	40056-T	Purchase 40056-T
CF®640R dCTP 25 nmol	CF®640R 40066	642/662 nm Purchase 40066	5 nmol	40066-T	Purchase 40066-T
CF®647 dCTP 25 nmol	CF®647 40028	650/665 nm Purchase 40028	5 nmol	40028-T	Purchase 40028-T
CF®660R dCTP 25 nmol	CF®660R 40068	663/682 nm Purchase 40068	5 nmol	40068-T	Purchase 40068-T

1. CF®594 dCTP is not compatible with PCR labeling of DNA probes.

CF is a registered trademark of Biotium, Inc. Alexa Fluor, Texas Red, and DyLight are registered trademarks of Thermo Fisher Scientific.

References

Download a list of curated [CF® Dye references](#).

This datasheet was generated on June 13, 2026 at 07:32:18 AM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/dctp-lyophilized-powder-cf-dye-conjugates/>