dUTP CF® Dye Conjugates

Labeled with our superior CF® dyes, dUTP conjugates can be used for TUNEL assay or to synthesize labeled DNA probes for in situ hybridization and nucleic acid blotting applications.



Product attributes

Call us: 800-304-5357

Colors

Blue, Green, Orange, Red, Far-red, Near-infrared

Email: btinfo@biotium.com

Product Description

CF® dye-dUTP can be used for TUNEL assay, microarray, or to synthesize labeled DNA probes for in situ hybridization and nucleic acid 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 8 CF® dve colors from blue to near infrared
- Biotium's CF® dyes have superior brightness, photostability, and are water-soluble

Supplied in lyophilized form. For PCR applications, Taq polymerase should be used with dUTP conjugates, because dUTP inhibits archaeal polymerases such as Pfu and Vent®. Also see our CF® Dye dCTP 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.

Note: Conjugates of blue-fluorescent dyes like CF®350, CF®405S and CF®405M are not recommended for detecting low abundance targets and may be challenging to use in tissue specimens. Blue dyes have lower fluorescence and photostability, and cells and tissue have high autofluorescence in blue wavelengths, resulting in lower signal to noise compared to other colors.

dUTP CF® Dye Conjugates

Product	Dye	Ex/Em	Size	SKU	Purchase
CF®405S dUTP ¹	CF®405S	404/431 nm	5 nmol	40004-T	Purchase 40004-T
			25 nmol	40004	Purchase 40004
CF®405M dUTP	CF®405M	408/452 nm	5 nmol	40100-T	Purchase 40100-T
			25 nmol	40100	Purchase 40100
CF®488A dUTP	CF®488A	490/515 nm	5 nmol	40008-T	Purchase 40008-T
			25 nmol	40008	Purchase 40008
CF®543 dUTP	CF®543	541/560 nm	5 nmol	40002-T	Purchase 40002-T
			25 nmol	40002	Purchase 40002
CF®568 dUTP	CF®568	562/583 nm	5 nmol	40005-T	Purchase 40005-T
			25 nmol	40005	Purchase 40005
CF®594 dUTP	CF®594	593/614 nm	5 nmol	40006-T	Purchase 40006-T
			25 nmol	40006	Purchase 40006
CF®640R dUTP	CF®640R	642/662 nm	5 nmol	40007-T	Purchase 40007-T
			25 nmol	40007	Purchase 40007
CF®680R dUTP ²	CF®680R	680/701 nm	5 nmol	40003-T	Purchase 40003-T
			25 nmol	40003	Purchase 40003

^{1.} CF®405S-dUTP may not be suitable for TUNEL staining in tissues due to blue autofluorescence in tissues and lower incorporation efficiency in tissue sections compared to other CF®dye dUTP conjugates. 2. CF®680R-dUTP has been tested in TUNEL staining of cells, but may not be efficiently incorporated in tissue sections.

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

References

Download a list of curated CF® Dye references.

This datasheet was generated on November 16, 2025 at 05:04:32 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/dutp-lyophilized-powder-cf-dye-conjugates/