

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

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

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

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 June 15, 2026 at 05:58:18 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/dutp-lyophilized-powder-cf-dye-conjugates/>