# **CF® Dye TCO**

CF® Dye with TCO (trans-cyclooctene) groups react with tetrazine-labeled molecules via copper-free click chemistry.



#### **Product attributes**

Chemical reactivity (reacts with)

Functional group

TCO (trans-cyclooctene)

Storage Conditions

Store at -10 to -35 °C, Protect from light

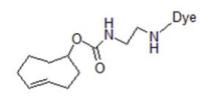
Call us: 800-304-5357 Email: btinfo@biotium.com

### **Product Description**

CF® Dye with TCO group reacts with tetrazine containing molecules via copper-free click chemistry. Copper-free bioorthogonal reactions are useful for labeling live cells, or when there are concerns about native protein function loss with copper in cell extracts. Membrane-permeant CF® Dye methyltetrazine options are also available for intracellular labeling reactions.

Available CF® Dye TCO derivatives: CF®488A, CF®532, CF®555, CF®568, CF®594, CF®640R, CF®647, CF®680, and CF®680R including membrane-permeant CF®500, and CF®650 TCO.

- Fluorescent labeling of tetrazine groups on target molecules via copper-free bioorthonal chemistry.
- Membrane-impermeant and membrane-permeant forms for surface or intracellular labeling of live cells or cell extracts.
- Bright, photostable and water-soluble CF® Dyes are excellent options for fluorescent labeling.



We also offer CF® Dye tetrazine and CF® Dye methyltetrazine.

#### 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. For more information download the CF® Dye Brochure.

## **CF Dye TCO: Product & SKU Summary**

Product	Conjugation	Ex/Em	Size	SKU	Purchase
CF®488A TCO	CF®488A	490/515 nm	1 mg	96040	Purchase 96040
CF®500 TCO	CF®500	500/510 nm	1 mg	96041	Purchase 96041
CF®532 TCO	CF®532	527/558 nm	1 mg	96042	Purchase 96042
CF®555 TCO	CF®555	555/565 nm	1 mg	96043	Purchase 96043
CF®568 TCO	CF®568	562/583 nm	1 mg	96044	Purchase 96044
CF®594 TCO	CF®594	593/614 nm	1 mg	96045	Purchase 96045
CF®640R TCO	CF®640R	642/662 nm	1 mg	96046	Purchase 96046
CF®647 TCO	CF®647	650/665 nm	1 mg	96047	Purchase 96047
CF®650 TCO	CF®650	650/670 nm	1 mg	96048	Purchase 96048
CF®680 TCO	CF®680	681/698 nm	1 mg	96051	Purchase 96051
CF®680R TCO	CF®680R	680/701 nm	1 mg	96049	Purchase 96049

This datasheet was generated on November 3, 2025 at 12:11:35 PM. Visit product page to check for updated information before use. Product link: <a href="https://biotium.com/product/cf-dye-tco/">https://biotium.com/product/cf-dye-tco/</a>