

## CF® Dye TCO

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



### Product attributes

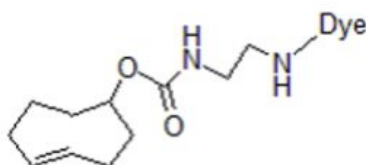
Chemical reactivity (reacts with)	Tetrazine
Functional group	TCO (trans-cyclooctene)
Storage Conditions	Store at -10 to -35 °C, Protect from light

## 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 bioorthogonal 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	<a href="#">Purchase 96040</a>
CF®500 TCO	CF®500	500/510 nm	1 mg	96041	<a href="#">Purchase 96041</a>
CF®532 TCO	CF®532	527/558 nm	1 mg	96042	<a href="#">Purchase 96042</a>
CF®555 TCO	CF®555	555/565 nm	1 mg	96043	<a href="#">Purchase 96043</a>
CF®568 TCO	CF®568	562/583 nm	1 mg	96044	<a href="#">Purchase 96044</a>
CF®594 TCO	CF®594	593/614 nm	1 mg	96045	<a href="#">Purchase 96045</a>
CF®640R TCO	CF®640R	642/662 nm	1 mg	96046	<a href="#">Purchase 96046</a>
CF®647 TCO	CF®647	650/665 nm	1 mg	96047	<a href="#">Purchase 96047</a>
CF®650 TCO	CF®650	650/670 nm	1 mg	96048	<a href="#">Purchase 96048</a>
CF®680 TCO	CF®680	681/698 nm	1 mg	96051	<a href="#">Purchase 96051</a>
CF®680R TCO	CF®680R	680/701 nm	1 mg	96049	<a href="#">Purchase 96049</a>

This datasheet was generated on December 20, 2025 at 06:52:17 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/cf-dye-tco/>