

CF® Dye Methyltetrazine

CF® Dye with methyltetrazine reacts with cyclooctene (TCO, methylcyclopropene) labeled molecules via copper-free click chemistry.



Product attributes

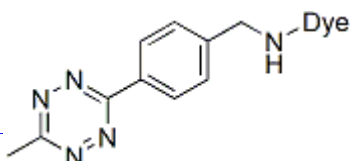
Chemical reactivity (reacts with)	Cyclooctene (TCO, methylcyclopropene) groups
Functional group	Tetrazine/Methyltetrazine
Storage Conditions	Store at -10 to -35 °C, Protect from light

Product Description

CF® Dye with a methyltetrazine group react with cyclooctene (TCO, methylcyclopropene) containing molecules via copper-free click chemistry. This copper-free bioorthogonal reaction is 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 methyltetrazine 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 methyltetrazine.

- Fluorescent labeling of cyclooctene (TCO, methylcyclopropene) 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.



CF® Dye methyltetrazines have improved chemical stability compared to [CF® Dye tetrazines](#). We also offer [CF® Dye TCO](#).

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 Methyltetrazine: Product & SKU Summary

Product	Conjugation	Ex/Em	Size	SKU	Purchase
CF@488A Methyltetrazine	CF@488A	490/515 nm	1 mg	96028	Purchase 96028
CF@500 Methyltetrazine	CF@500	500/510 nm	1 mg	96029	Purchase 96029
CF@532 Methyltetrazine	CF@532	527/558 nm	1 mg	96030	Purchase 96030
CF@555 Methyltetrazine	CF@555	555/565 nm	1 mg	96031	Purchase 96031
CF@568 Methyltetrazine	CF@568	562/583 nm	1 mg	96032	Purchase 96032
CF@594 Methyltetrazine	CF@594	593/614 nm	1 mg	96033	Purchase 96033
CF@640R Methyltetrazine	CF@640R	642/662 nm	1 mg	96034	Purchase 96034
CF@647 Methyltetrazine	CF@647	650/665 nm	1 mg	96035	Purchase 96035
CF@650 Methyltetrazine	CF@650	650/670 nm	1 mg	96036	Purchase 96036
CF@680 Methyltetrazine	CF@680	681/698 nm	1 mg	96039	Purchase 96039
CF@680R Methyltetrazine	CF@680R	680/701 nm	1 mg	96037	Purchase 96037