# **CF® Dye Methyltetrazine**

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



### Product attributes

Call us: 800-304-5357

Chemical reactivity (reacts with)

Functional group

Tetrazine/Methyltetrazine

Storage Conditions

Cyclooctene (TCO, methylcyclopropene) groups

Tetrazine/Methyltetrazine

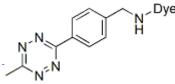
Email: btinfo@biotium.com

### **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 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.



CF® Dye methyltetrazines have improved chemical stability compared to <u>CF® Dye tetrazines</u>. We also offer <u>CF® Dye TCO</u>.

#### 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 |

This datasheet was generated on November 3, 2025 at 12:11:07 PM. Visit product page to check for updated information before use. Product link: <a href="https://biotium.com/product/cf-dye-methyltetrazine/">https://biotium.com/product/cf-dye-methyltetrazine/</a>