

## CF® Dye Tetrazine

CF® Dyes with a tetrazine group react with cyclooctene (TCO, methylcyclopropene) labeled molecules via copper-free click chemistry.



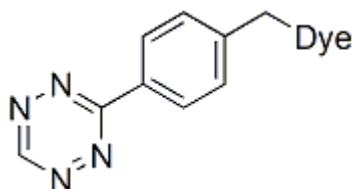
### Product attributes

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

## Product Description

Tetrazine CF® Dyes react with cyclooctene (TCO, methylcyclopropene) labeled molecules via copper-free click chemistry. These copper-free bioorthogonal reactions are useful for labeling fixed cells or when there are concerns about native protein function loss with copper in cell extracts.

- Fluorescent labeling of cyclooctene (TCO, methylcyclopropene) groups on target molecules via copper-free bioorthogonal chemistry.
- Faster reaction kinetics than methyltetrazine derivatives.
- Bright, photostable, and water-soluble CF® Dyes are excellent options for fluorescent labeling.



We also offer [CF® Dye Methyltetrazines](#) and [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 Tetrazine

Dye	Ex/Em	Size	Catalog No.	Dye Features
<a href="#">CF@488A</a>	490/516 nm	1 mg	<a href="#">96054</a>	<a href="#">CF@488A Features</a>
<a href="#">CF@568</a>	562/584 nm	1 mg	<a href="#">96055</a>	<a href="#">CF@568 Features</a>
<a href="#">CF@647</a>	650/668 nm	1 mg	<a href="#">96056</a>	<a href="#">CF@647 Features</a>
<a href="#">CF@660C</a>	667/685 nm	1 mg	<a href="#">96096</a>	<a href="#">CF@660C Features</a>
<a href="#">CF@680</a>	681/698 nm	1 mg	<a href="#">96097</a>	<a href="#">CF@680 Features</a>

Alexa Fluor and DyLight are registered trademarks of Thermo Fisher Scientific.

This datasheet was generated on January 9, 2026 at 12:57:30 AM. Visit product page to check for updated information before use.  
Product link: <https://biotium.com/product/cf-dye-tetrazine/>