

Revised: March 14, 2019

## **Product Information**

# **CF® Dye Methyltetrazine or Biotin Methylterazine**

Unit Size: 1.0 mg

#### **Technical Summary**

Cat. No.	CF® Dye	Abs <sub>max</sub> (nm)	Em <sub>max</sub> (nm)	Extinction coefficient	MW
96028	CF®488A	490	515	70,000	~1097
96029	CF®500	500	510	70,000	~475
96030	CF®532	527	558	96,000	~869
96031	CF®555	555	565	150,000	~1142
96032	CF®568	562	583	100,000	~897
96033	CF®594	593	614	115,000	~912
96034	CF®640R	642	662	105,000	~1015
96035	CF®647	650	665	240,000	~1168
96036	CF®650	650	670	100,000	~657
96037	CF®680R	680	701	140,000	~1095
96038	Biotin				~427
96039	CF®680	681	698	210,000	~3336

### Storage and Handling

Store CF® dye methyltetrazine or biotin methylterazine at -20°C, protected from light. Product is stable for at least 12 months from date of receipt if stored as recommended. Stock solution in DMSO can be stored at  $\leq$  -20°C for at least 12 months

#### **Product Description**

For this set of products, we provide our CF® dyes or biotin in the methyltetrazine form. CF® dyes are exceptionally bright and photostable, making it the perfect dye for fluorescence detection. CF® dye with methyltetrazine reacts with cyclooctene (TCO, methylcyclopropene) labeled molecule via copper free click chemistry. This copper free bioorthogonal reaction allows staining the surface of live cells or having concerns about native protein function loss with copper in cell extracts.

CF®500 methyltetrazine and CF®650 methyltetrazine are designed to use for intracellular copper free reaction with cyclooctene.

#### Other Related Products

You may also be interested in the following related products from Biotium:

- · A full selection CF® reactive dyes and CF® dye conjugates.
- CF® dye-labeled azides.
- CF® dye-labeled alkyne.
- CF® dye streptavidin.
- · CF® dye BCN.

Please visit www.biotium.com to view our full selection of innovative products for life science research.

CF® dye technology is covered by pending U.S. and international patents. Materials from Biotium are sold for research use only, and are not intended for food, drug, household, or cosmetic use.