



# **Product Information**

## **CF™ Dye TCO or Biotin TCO**

Unit Size: 1.0 mg

### **Technical Summary**

Cat. No.	CF™ Dye	Abs <sub>max</sub> (nm)	Em <sub>max</sub> (nm)	Extinction coefficient	MW
96040	CF™488A	490	515	70,000	~1108
96041	CF™500	500	510	70,000	~486
96042	CF™532	527	558	96,000	~880
96043	CF™555	555	565	150,000	~1153
96044	CF™568	562	583	100,000	~908
96045	CF™594	593	614	115,000	~923
96046	CF™640R	642	662	105,000	~1026
96047	CF™647	650	665	240,000	~1179
96048	CF™650	650	670	100,000	~668
96049	CF™680R	680	701	140,000	~1106
96050	Biotin				~438
96051	CF™680	681	698	210,000	~3347

#### Storage and Handling

Store  $CF^{TM}$  dye TCO or biotin TCO 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**

CF™ dye or biotin with TCO (trans-cyclooctene) reacts with tetrazine 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™ dyes are exceptionally bright and photostable, making it the perfect dye for fluorescence detection.

CF™500 TCO and CF™650 TCO are designed to use for intracellular copper free reaction with tetrazine.

#### **Other Related Products**

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

- A full selection CF<sup>™</sup> reactive dyes and CF<sup>™</sup> dye conjugates
- CF<sup>™</sup> dye-labeled azides.
- CF<sup>™</sup> dye-labeled alkyne.
- CF™ dye streptavidin.
- CF™ dye methylterazine.
- 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.