

Product Information

CF™ Dye Tetrazine

Unit Size: 1.0 mg

Technical Summary

Cat. No.	CF™ Dye	Abs _{max} (nm)	Em _{max} (nm)	Extinction coefficient	MW
96054	CF™488A	490	515	70,000	~1083
96055	CF™568	562	583	100,000	~883
96056	CF™647	650	665	240,000	~1154

Storage and Handling

Store CF™ dye tetrazine 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 ≤ -20°C for at least 12 months.

Product Description

For this set of products, we provide our CF™ dyes in the tetrazine form. CF dyes are exceptionally bright and photostable, making it the perfect dye for fluorescence detection. CF™ dye with tetrazine 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™ dyes tetrazine allow faster reaction kinetics than CF™ dyes methyltetrazine.

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.