

May 31, 2013

Product Information

MTS-CF™ Dye

Unit Size: 1 mg

Technical Summary

Cat. No.	CF™ Dye	Abs _{max} (nm)	Em _{max} (nm)	Extinction coefficient	MW
92096	CF™640R	642	662	105,000	~969
92097	CF™488A	490	515	70,000	~1051
92098	CF™543	541	560	100,000	~1024
92099	CF™594	593	614	115,000	~866

Storage and Handling

Store MTS-CF□ dye at -20°C, protected from light. Product is stable for at least 12 months from date of receipt if stored as recommended. Stock solutions may be prepared in DMSO and can be stored at ≤ -20°C for at least 12 months.

Product Description

MTS-CF™ dye is a fast-reacting and highly selective thiol-reactive label that permits one to attach CF™ dye to a thiol group. CF™ dyes are bright, photostable and water-soluble, making them an excellent choice for fluorescent labeling.

Related Products

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

- CF™ dye succinimidyl (NHS) ester, hydrazide, maleimide, aminooxy, alkyne, and azide derivatives
- CF™ antibody conjugates and protein labeling kits
- . Mix-n-stain CF™ antibody labeling kits

Please visit our website at www.biotium.com for details.

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.