

CF® Dye MTS

CF® dye MTS are fast-reacting, highly selective thiol-reactive derivatives for reversible labeling of thiol groups with fluorescent CF® dyes.



Product attributes

Chemical reactivity (reacts with)	Thiols
Functional group	MTS
Storage Conditions	Store at -10 to -35 °C or below, Protect from light

Product Description

CF® dyes with a methanethiosulfonate (MTS) group are fast-reacting and highly selective thiol-reactive derivatives that can be used to attach fluorescent CF® dyes to thiol groups. Labeling can be easily reversed using reducing agents like DTT or TCEP.

- Rapid and highly-selective labeling of thiol groups on target molecules with fluorescent CF® dyes.
- Labeling is reversible using reducing reagents like TCEP and DTT.
- Bright, photostable and water-soluble CF® dyes are excellent options for fluorescent labeling.

Superior CF® Dyes

CF® dyes are Biotium's line of next-generation fluorescent dyes that have improved brightness, photostability and water solubility compared to other commercially available fluorescent dyes.

Learn more about [CF® Dyes](#). For more information download the [CF® Dye Brochure](#).

Product	Conjugation	Ex/Em	Size	Catalog No.	Purchase
CF®488A MTS	CF®488A	490/515 nm	1 mg	92097	Purchase 92097
CF®543 MTS	CF®543	541/560 nm	1 mg	92098	Purchase 92098
CF®594 MTS	CF®594	593/614 nm	1 mg	92099	Purchase 92099
CF®640R MTS	CF®640R	642/662 nm	1 mg	92096	Purchase 92096

References

Download a list of [CF® dye references](#).

This datasheet was generated on June 16, 2026 at 10:19:51 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/cf-dye-mts/>