

Product Information

CF™ Dye Amine

Unit Size: 1 mg

Technical Summary

Cat. No.	CF™ Dye	Abs _{max} (nm)	Em _{max} (nm)	Extinction coefficient	MW
92035	CF™350	347	448	18,000	~537
92036	CF™405S	408	452	33,000	~725
92037	CF™488A	490	515	70,000	~597
92038	CF™555	555	565	150,000	~943
92039	CF™568	562	583	100,000	~756
92040	CF™594	593	614	115,000	~771
92041	CF™633	630	650	100,000	~863
92043	CF™640R	642	662	105,000	~1034
92042	CF™647	650	665	240,000	~969
96010	CF™660R	663	682	100,000	~930
92102	CF™750	755	777	250,000	~2963
92065	CF™770	770	797	250,000	~3180

Storage and Handling

Store CF™ dye amine 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 or dH₂O and can be stored at ≤ -20°C for at least 12 months.

Product Description

CF™ dye amines can be conjugated to activated carboxylic acids in proteins or other molecules. Our CF™ amine derivatives are bright 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, aminoxy, alkyne, and azide derivatives
- CF™ antibody conjugates and protein 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.