

## Product Information

### CF™568, Hydrazide

Catalog Number **SCJ4600025**

Storage Temperature  $-20\text{ }^{\circ}\text{C}$

## TECHNICAL BULLETIN

### Product Description

CF™568, hydrazide can be used as a fixable polar tracer for visualizing cell morphology, or for labeling biomolecules with an aldehyde or ketone group (such as carbohydrate molecules after peroxidation with periodate).

CF568 is a red fluorescent dye with an excitation spectrum optimally matching the 568 nm line of the Ar-Kr mixed-gas laser. Antibody conjugates of CF568 are much brighter than those of Alexa Fluor® 568. In addition, the photostability of CF568 is superior to that of Alexa Fluor 568, making CF568 a much better choice for demanding applications such as confocal microscopy and single molecule imaging.

#### CF568 dye properties:

Abs/Em Maxima: 562/583 nm (See Figure 1)

Extinction coefficient: 100,000

Molecular weight: ~700

$A_{280}/A_{\text{max}}$  or CF (correction factor for estimating degree of protein labeling): 0.08

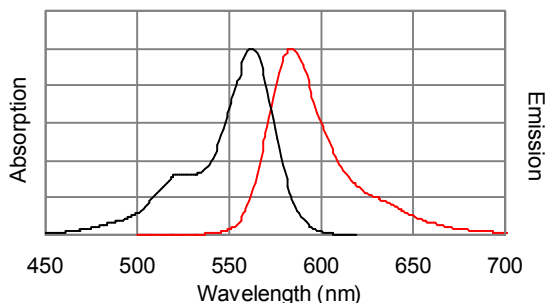
Flow cytometry laser line: 488, 532, or 561 nm

Microscopy laser line: 568 nm

Direct replacement for: Alexa Fluor 568 and Rhodamine Red

### Figure 1.

Absorption and emission spectra of CF568 conjugated to goat anti-mouse IgG in PBS.



### Precautions and Disclaimer

This product is for R&D use only, not for drug, household, or other uses. Please consult the Material Safety Data Sheet for information regarding hazards and safe handling practices.

### Preparation Instructions

Stock solutions may be prepared in water or PBS.

Stock solutions may be stored at  $-20\text{ }^{\circ}\text{C}$  for at least 6 months.

### Storage/Stability

Store the dye desiccated at  $-20\text{ }^{\circ}\text{C}$ . When stored as directed, the dye should remain active for at least 6 months.

This product is distributed by Sigma-Aldrich Co. under the authorization of Biotium, Inc. This product is covered by one or more US patents and corresponding patent claims outside the US patents or pending applications owned or licensed by Biotium, Inc. including without limitation: 12/334,387; 12/607,915; 12/699,778; 12/850,578; 61/454,484. In consideration of the purchase price paid by the buyer, the buyer is hereby granted a limited, non-exclusive, non-transferable license to use only the purchased amount of the product solely for the buyer's own internal research in a manner consistent with the accompanying product literature. Except as expressly granted herein, the sale of this product does not grant to or convey upon the buyer any license, expressly, by implication or estoppel, under any patent right or other intellectual property right of Biotium, Inc.

Buyer shall not resell or transfer this product to any third party, or use the product for any commercial purposes, including without limitation, any diagnostic, therapeutic or prophylactic uses. This product is for research use only. Any other uses, including diagnostic uses, require a separate license from Biotium, Inc. For information on purchasing a license to use this product for purposes other than research, contact Biotium, Inc., 3159 Corporate Place, Hayward, CA 94545, Tel: (510) 265-1027. Fax: (510) 265-1352. Email: [btinfo@biotium.com](mailto:btinfo@biotium.com).

CF is a trademark of Biotium.

Alexa Fluor is a registered trademark of Invitrogen;

AKN,MAM 10/11-1