

Product Information

CF™680, Free Acid

Catalog Number: 92070

Unit Size: 1 mg

Color and Form: Dark blue solid.

Storage and Handling

Store CF™680 free acid desiccated at $\leq -20^{\circ}\text{C}$. When stored as directed, CF™680 free acid should be stable for at least 12 months from the time of receipt.

Technical Summary

Abs/Em Maxima: 681/698 nm (See Figure 1)

Extinction coefficient: 210,000

Molecular weight: ~3241

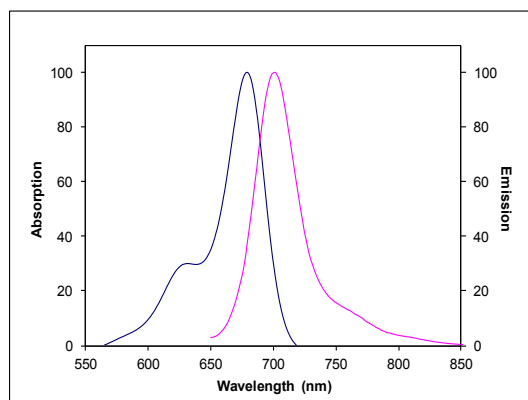


Figure 1. Absorption and emission spectra of CF™680 Free Acid in water

Product Application

CF™680 free acid has no reactive group and cannot be used for labeling. CF680 free acid could be used for the following applications:

In vitro cell-based assays

As a control for cell-based assays that monitor binding of a dye-labeled agent

- Validation of optical agent for in vivo administration
- Evaluation of binding specificity

In vivo imaging

For evaluation of behavior and clearance of the dye itself

- Timing of dye clearance from the animal's body
- Retention of dye in certain organs or sites (e.g., liver or kidneys)

Other Related Products

Cat. #	Product Name	Unit Size
92139	CF680 succinimidyl ester	1 umol
92240	Mix-n-Stain CF680 antibody labeling kit, 1X(50-100ug) labeling	1 labeling
92262	Mix-n-Stain CF680 antibody labeling kit, 1X(20-50ug) labeling	1 labeling
92282	Mix-n-Stain CF680 antibody labeling kit, 1X(5-20ug) labeling	1 labeling
92160	VivoBrite rapid antibody labeling kit for small animal in vivo imaging, CF680 SE	3 labelings
29007	Annexin V, CF680 conjugate, lyophilized	25 ug
92220	CF680 SE protein labeling kit	3 labelings
20063-1	CF680 Goat anti-mouse IgG (H+L), F(ab') ₂ fragment, 2 mg/mL	50 uL
20065-1	CF680 Goat anti-mouse IgG (H+L), highly cross-adsorbed, 2 mg/mL	50uL
20064-1	CF680 Goat anti-rabbit IgG (H+L), F(ab') ₂ fragment, 2 mg/mL	50uL
20067-1	CF680 Goat anti-rabbit IgG (H+L), highly cross-adsorbed, 2 mg/mL	50uL