

Product Information

CF™770, Free Acid

Catalog Number: 92066

Unit Size: 1 mg

Color and Form: Dark blue solid.

Storage and Handling

Store CF™770 free acid desiccated at $\leq -20^{\circ}\text{C}$. When stored as directed, CF™770 free acid should be stable for at least 12 months from the time of receipt.

Technical Summary

Abs/Em Maxima: 770/797 nm (See Figure 1)

Extinction coefficient: 220,000

Molecular weight: ~3092

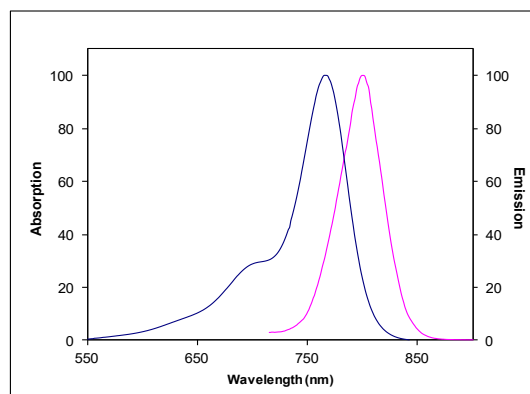


Figure 1. Absorption and emission spectra of CF™770 Free Acid in water

Product Application

CF™770 free acid has no reactive group and cannot be used for labeling. CF770 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
92150	CF770 succinimidyl ester	1 umol
92242	Mix-n-Stain CF770 antibody labeling kit, 1X(50-100ug) labeling	1 labeling
92265	Mix-n-Stain CF770 antibody labeling kit, 1X(20-50ug) labeling	1 labeling
92285	Mix-n-Stain CF770 antibody labeling kit, 1X(5-20ug) labeling	1 labeling
92162	VivoBrite rapid antibody labeling kit for small animal in vivo imaging, CF770 SE	3 labelings
29046	Annexin V, CF770 conjugate, lyophilized	25 ug
92222	CF770 SE protein labeling kit	3 labelings
20077-1	CF770 Goat Anti-Mouse IgG (H+L) (min X Human, Bovine, Horse, Rabbit, and Swine), 2 mg/mL	50 uL
20078-1	CF770 Goat Anti-Rabbit IgG (H+L) (min X Human, Mouse, and Rat), 2 mg/mL	50uL