

Product Information

CF™790, Free Acid

Catalog Number: 92159

Unit Size: 0.25 μ mole

Color and Form: Dark blue solid.

Storage and Handling

Store CF™790 free acid desiccated at $\leq -20^{\circ}\text{C}$. When stored as directed, CF™790 free acid should be stable for at least 12 months from the time of receipt.

Technical Summary

Abs/Em Maxima: 784/806 nm (See Figure 1)

Extinction coefficient: 210,000

Molecular weight: ~3221

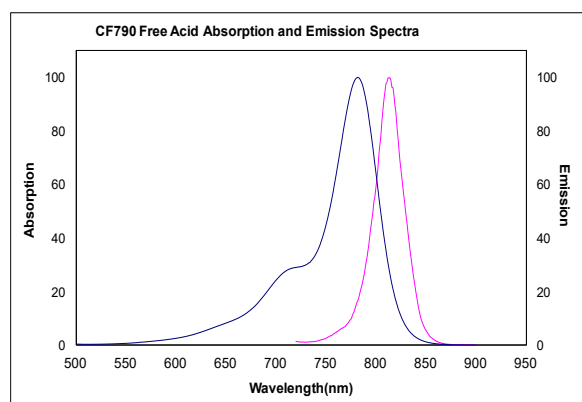


Figure 1. Absorption and emission spectra of CF™790 Free Acid in water

Product Application

CF™790 free acid has no reactive group and cannot be used for labeling. CF790 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
29047	Annexin V, CF790 conjugate, lyophilized	25 μ g
20345-50 μ L	CF790 Donkey Anti-Goat IgG (H+L), highly cross-adsorbed, 2mg/mL	50 μ L
20363-50 μ L	CF790 Donkey Anti-Mouse IgG (H+L), highly cross-adsorbed, 2mg/mL	50 μ L
20344-50 μ L	CF790 Donkey Anti-Rabbit IgG (H+L), highly cross-adsorbed, 2mg/mL	50 μ L
92155	CF790, succinimidyl ester	0.25 μ mol
20378-50 μ L	CF790 Goat Anti-Mouse IgG (H+L), 2 mg/mL	50 μ L
20342-50 μ L	CF790 Goat Anti-Mouse IgG (H+L), highly cross-adsorbed, 2 mg/mL	50 μ L
20379-50 μ L	CF790 Goat Anti-Rabbit IgG (H+L), 2 mg/mL	50 μ L
20343-50 μ L	CF790 Goat Anti-Rabbit IgG (H+L), highly cross-adsorbed, 2 mg/mL	50 μ L
92288	CF790 Mix-n-Stain Antibody Labeling Kit	1 X 5-20 μ g labeling
92163	VivoBrite rapid antibody labeling kit for small animal in vivo imaging, CF790 SE	3 labelings