

# Product Information

## CF™750, Free Acid

**Catalog Number:** 92071

**Unit Size:** 1 mg

**Color and Form:** Dark blue solid.

### Storage and Handling

Store CF™750 free acid desiccated at ≤ -20°C. When stored as directed, CF™750 free acid should be stable for at least 12 months from the time of receipt.

### Technical Summary

**Abs/Em Maxima:** 755/777 nm (See Figure 1)

**Extinction coefficient:** 250,000

**Molecular weight:** ~2963

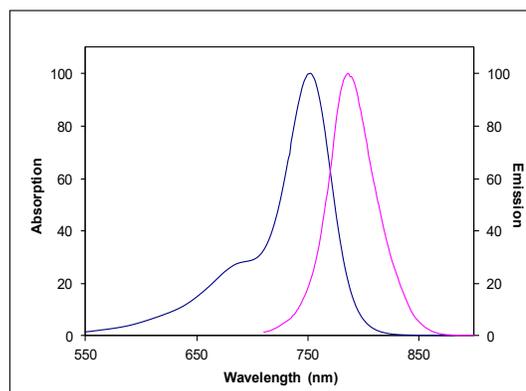


Figure 1. Absorption and emission spectra of CF™750 Free Acid in water

### Product Application

CF™750 free acid has no reactive group and cannot be used for labeling. CF750 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
92142	CF750 succinimidyl ester	1 umol
92241	Mix-n-Stain CF750 antibody labeling kit, 1X(50-100ug) labeling	1 labeling
92264	Mix-n-Stain CF750 antibody labeling kit, 1X(20-50ug) labeling	1 labeling
92284	Mix-n-Stain CF750 antibody labeling kit, 1X(5-20ug) labeling	1 labeling
92161	VivoBrite rapid antibody labeling kit for small animal in vivo imaging, CF750 SE	3 labelings
29006	Annexin V, CF750 conjugate, lyophilized	25 ug
92221	CF750 SE protein labeling kit	3 labelings
20070-1	CF750 Goat Anti-Mouse IgG (H+L), 2 mg/mL	50 uL
20073-1	CF750 Goat Anti-Rabbit IgG (H+L), 2 mg/mL	50uL
20362-50uL	CF750 Donkey Anti-Goat IgG (H+L), highly cross-adsorbed, 2mg/mL	50uL
20298-1	CF750 Donkey Anti-Rabbit IgG (H+L), highly cross-adsorbed, 2mg/mL	50uL