

Product Information

CF Dye-Dextran 250,000 MW, Anionic & Fixable

Catalog Number:

| Cat. No. | Conjugate | Abs _{max} (nm) | Em _{max} (nm) |
|----------|-----------|-------------------------|------------------------|
| 80134 | CF™488A | 490 | 515 |
| 80135 | CF™680 | 681 | 698 |
| 80136 | CF™750 | 755 | 777 |
| 80125 | CF™770 | 770 | 797 |

Unit Size: 1 mg

Storage and Handling

Store at -20 °C and protect from light, especially when in solution. Product is stable for at least 1 year from date of receipt when stored as recommended.

Spectral Properties

See table above.

Preparation of Stock Solution

Dextran conjugates are generally soluble in water and aqueous buffers to at least 10 mg/mL. Aqueous solutions of dextrans can be stored at 2-6 °C for several weeks, with the addition of sodium azide to a final concentration of 2 mM to inhibit bacterial growth. For long-term or preservative-free storage, divide the aqueous solution into aliquots and freeze at -20 °C.

Product Description

CF dye labeled dextran 250,000 MW could be used as a fluorescent probe to study cellular processes such as cell permeability, phagocytosis, endocytosis, or to study mechanisms of biomolecular delivery. CF™ dyes are superior fluorescent dyes with exceptional brightness and remarkable photostability.

Related Products

You may also be interested in the following related products from Biotium:

- CF™ dye succinimidyl (NHS) ester, hydrazide, maleimide, aminoxy, alkyne, and azide derivatives
- CF™ dye conjugates including antibodies, streptavidin, phalloidin, transferrin, cholera toxin, BSA, lectins, nucleotides, and more
- Antibody and protein labeling kits

Please visit our website at www.biotium.com for details.

CF dye technology is covered by pending U.S. and international patents. Materials from Biotium are sold for research use only, and are not intended for food, drug, household, or cosmetic use.