

www.biotium.com September 28, 2017

# **Product Information**

## CF Dye-Dextran 3,500 MW, Anionic & **Fixable**

#### **Catalog Number:**

Cat. No.	Conjugate	Abs <sub>max</sub> (nm)	Em <sub>max</sub> (nm)
80137	CF™350	347	448

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 3,500 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, aminooxy, alkyne, and azide derivatives
- CF™ dye conjugates including antibodies, streptavidin, phalloidin, transferrin, cholera toxin, BSA, lectins, nucleotides, and more
- · Antibody and protein labeling kits
- CF dye dextrans, anionic and fixable (10,000MW, 40,000MW, 70,000MW, 150,000MW and 250,000MW).

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.