

wing Products for Science™ www.biotium.com

Revised: November 19. 2021

Product Information

CF® Dye Dextran 70,000 MW, Anionic & Fixable

Product List

Catalog no.	CF® Dye	Ex/Em (nm)
80138	CF®350	347/448
80117	CF®488A	490/515
80141	CF®633	630/650
80129	CF®680	681/698
80130	CF®750	755/777
80123	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 70,000 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

Catalog number	Product
80137	CF® Dye Dextran 3,500 MW, Anionic and Fixable
80110 - 80121	CF® Dye Dextran 10,000 MW, Anionic and Fixable
80122 - 80128	CF® Dye Dextran 40,000 MW, Anionic and Fixable
80124 - 80133	CF® Dye Dextran 150,000 MW, Anionic and Fixable
80125 - 80136	CF® Dye Dextran 250,000 MW, Anionic and Fixable
92004- 92007	CF® Dye Biocytin
92151- 96025	CF® Dye Hydrazide
00068- 00079	Cholera Toxin Subunit B CF® Dye Conjugates
00081- 00087	Transferrin (Human) CF® Dye Conjugates
00002 00026	α-Bungarotoxin, CF® Dye and Other Conjugates
20289- 20292	CF® Dye Bovine Serum Albumin (BSA)
92230- 92339	Mix-n-Stain™ CF® Dye Antibody Labeling Kits
92454 - 92411	Mix-n-Stain™ Maxi Antibody Labeling Kits, 1 mg Labeling
30068 - 30050	ViaFluor® SE Cell Proliferation Kits

Please visit our website at www.biotium.com for information on our life science research products, including environmentally friendly EvaGreen® qPCR master mixes, fluorescent CF® Dye antibody conjugates and reactive dyes, apoptosis reagents, fluorescent probes, and kits for cell biology research.

CF® Dye technology is covered by 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.