

CF® Dye Dextran 10,000 MW, Anionic and Fixable



CF® Dye labeled dextrans could be used as a fluorescent fluid-phase markers to study cell permeability, endocytosis, or mechanisms of biomolecular delivery. The dextran is 10,000 MW, and contains a formaldehyde-fixable free-amine group.

Product attributes

Probe cellular localization	Fluid phase tracer
Cell permeability	Membrane impermeant
Colors	Green, Orange, Red, Far-red, Near-infrared

Product Description

CF® Dye labeled dextrans could be used as a fluorescent fluid-phase markers to study cell permeability, endocytosis, or mechanisms of biomolecular delivery. The dextran is 10,000 MW, and contains a formaldehyde-fixable free-amine group.

CF® Dye Dextran 10,000 MW

Conjugation	Ex/Em	Size	Catalog No.	Dye Features
CF@488A	490/515 nm	1 mg	80110	CF@488A Features
CF@543	541/560 nm	1 mg	80111	CF@543 Features
CF@555	555/565 nm	1 mg	80112	CF@555 Features
CF@568	562/583 nm	1 mg	80113	CF@568 Features
CF@594	593/614 nm	1 mg	80114	CF@594 Features
CF@640R	642/662 nm	1 mg	80115	CF@640R Features
CF@680	681/698 nm	1 mg	80118	CF@680 Features
CF@680R	680/701 nm	1 mg	80116	CF@680R Features
CF@750	755/777 nm	1 mg	80119	CF@750 Features
CF@770	770/797 nm	1 mg	80120	CF@770 Features
CF@790	784/806 nm	1 mg	80121	CF@790 Features

References

Download a list of [CF® dye references](#).

This datasheet was generated on June 4, 2026 at 10:40:14 AM. Visit product page to check for updated information before use.
 Product link: <https://biotium.com/product/cf-dye-dextran-10000-mw-anionic-fixable/>