



Glowing products for science

Call us : [800-304-5357](tel:800-304-5357)

CF® Dye Biocytin

CF® Dye biocytin is a cell-impermeant, fixable polar tracer that is a conjugate of CF® Dye and biotin with an aldehyde-fixable primary amine. Polar tracers are commonly used to investigate cell-cell and cell-liposome fusion as well as membrane permeability and transport through gap junctions or cell uptake during pinocytosis.



Product attributes

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

Product Description

CF® Dye biocytin is a cell-impermeant, fixable polar tracer that is a conjugate of CF® Dye and biotin with an aldehyde-fixable primary amine. Polar tracers are commonly used to investigate cell-cell and cell-liposome fusion as well as membrane permeability and transport through gap junctions or cell uptake during pinocytosis.

CF® Dye biocytin is available in the following colors: CF@488A, CF@568, CF@594, and CF@640R.

CF® Dye Biocytin

CF® Dye Biocytin	Ex/Em	Size	Catalog No.	Dye Features
CF@488A	490/515 nm	0.5 mg	92004	CF@488A Features
CF@568	562/583 nm	0.5 mg	92005	CF@568 Features
CF@594	593/614 nm	0.5 mg	92006	CF@594 Features
CF@640R	642/662 nm	0.5 mg	92007	CF@640R Features

This datasheet was generated on June 15, 2026 at 01:06:15 AM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/cf-dye-biocytin/>