## **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**

Call us: 800-304-5357

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

Email: btinfo@biotium.com

## **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	<u>92004</u>	CF®488A Features
<u>CF®568</u>	562/583 nm	0.5 mg	<u>92005</u>	CF®568 Features
CF®594	593/614 nm	0.5 mg	<u>92006</u>	CF®594 Features
<u>CF®640R</u>	642/662 nm	0.5 mg	<u>92007</u>	CF®640R Features

This datasheet was generated on November 14, 2025 at 03:16:45 AM. Visit product page to check for updated information before use. Product link: <a href="https://biotium.com/product/cf-dye-bioctyin/">https://biotium.com/product/cf-dye-bioctyin/</a>