Glowing products for science

TMR Biocytin

TMR biocytin is a cell-impermeant, fixable polar tracer that combines the tetramethylrhodamine (TAMRA) with biotin and 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	Red
Excitation/Emission	544/571 nm
Conjugation	Tetramethylrhodamine (TRITC), Biotin

Product Description

TMR biocytin is a cell-impermeant, fixable polar tracer that combines the tetramethylrhodamine (TAMRA) with biotin and 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. TMR biocytin has proved to be an effective neuronal tracer in live tissue when applied by electroporation¹.

- λ_{Ex}/λ_{Em}: 544/571 nm
- Dark red solid soluble in DMF or DMSO
- MW: 871

We also offer CF® Dye-Biocytin, with a selection of our superior CF® dyes.

References

1. Journal of Neuroscience Methods (2015), http://dx.doi.org/doi:10.1016/j.jneumeth.2015.06.005

This datasheet was generated on November 20, 2025 at 08:03:04 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/tmr-biocytin/