

## 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

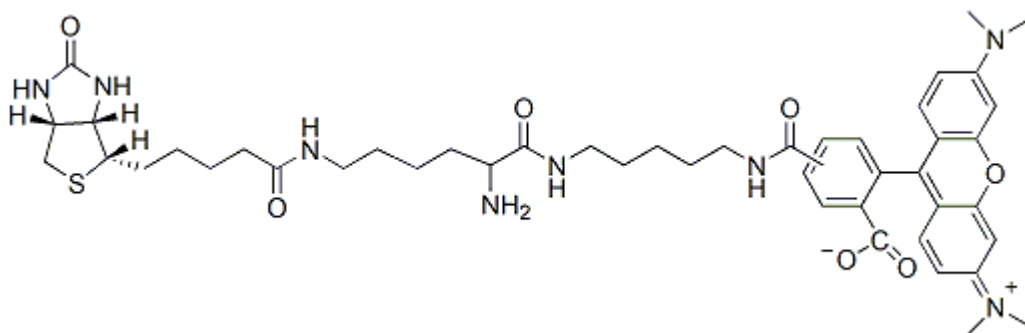
Probe cellular localization	Fluid phase tracer
Cell permeability	Membrane impermeant
Colors	Red
Excitation/Emission	544/571 nm
Conjugation	Tetramethylrhodamine (TRITC), Biotin

## 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 electroporation<sup>1</sup>.

- $\lambda_{Ex}/\lambda_{Em}$ : 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 January 18, 2026 at 10:51:28 AM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/tmr-biocytin/>