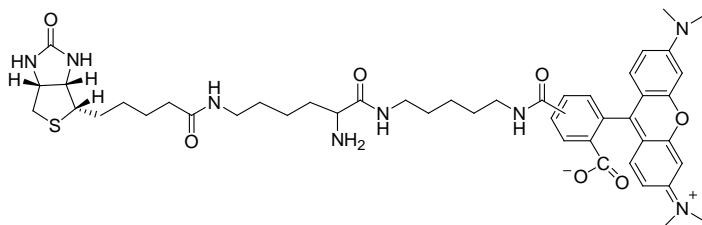


PRODUCT AND SAFETY DATA SHEET

PRODUCT NAME: TMR Biocytin**CATALOG #** 92003**MOLECULAR INFORMATION:** C₄₆H₈₂N₈O₇S
MWt: 871**PROPERTIES:****Color & Form**

Dark Red solid

Solubility

Soluble in DMF or DMSO

Absorption / Emission

544 / 571 nm

STORAGE AND HANDLING:

Store at -20°C

APPLICATION:

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 proved to be an effective neuronal tracer in live tissue when applied by electroporation¹.

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

TOXICITY:

Not established. Not listed by NTP, IARC or OSHA.

FIRST AID:

Potentially harmful. Avoid prolonged or repeated exposure. Avoid getting in eyes, on skin, or on clothing. Wash thoroughly after handling. If eye or skin contact occurs, wash affected areas with plenty of water for 15 minutes and seek medical advice. In case of inhaling or swallowing, move individual to fresh air and seek medical advice immediately.

Disclaimer: Materials from Biotium are sold for research use only, and are not intended for food, drug, household, or cosmetic use. Biotium is not liable for any damage resulting from handling or contact with this product.