

D-Biocytn TMR

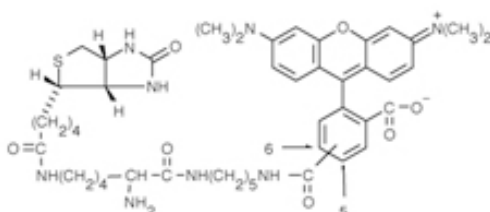
Catalog Number	Packaging Size
B011	5 mg

Storage upon receipt: -20°C

Introduction

The water-soluble neuronal tracer **5-(and-6)-tetramethylrhodamine biocytn (biocytn TMR)** can be detected in fixed tissues by using avidin or streptavidin probes. However, tetramethylrhodamine biocytn can also be observed directly in fixed or unfixed samples because the attached tetramethylrhodamine dye produces bright, orange-fluorescence (absorption/emission maxima ~555/580 nm).

Specifications

Molecular Formula:	C ₄₆ H ₆₀ N ₈ O ₇ S	
Molecular Weight:	869.09	
CAS Number:	-	
Storage Conditions:	-20°C	
Shipping Condition:	Room Temperature	

Applications

Water-soluble neuronal tracers

References:

[Evaluation of biotin-dye conjugates for use in an HPLC assay to assess relative binding of biotin derivatives with avidin and streptavidin.](#)

Wilbur DS, Pathare PM, Hamlin DK, Frownfelter MB, Kegley BB, Leung WY, Gee KR
Bioconjug Chem (2000) 11:584-598