

D-Biotin-4-fluorescein

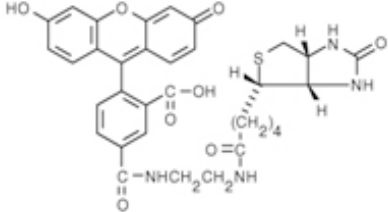
Catalog Number	Packaging Size
B009	10 mg

Storage upon receipt: -20°C

Introduction

Biotin-4-fluorescein can be used to quantitate available biotin-binding sites or the concentration of avidin or streptavidin through the strong quenching associated when this molecule binds to avidin.

Specifications

Molecular Formula:	C ₃₃ H ₃₂ N ₄ O ₈ S	
Molecular Weight:	644.70	
CAS Number:	-	
Storage Conditions:	-20°C	
Shipping Condition:	Room Temperature	

Applications

Quantitation of biotin-binding sites

References:

1. Evaluation of fluorescent compound interference in 4 fluorescence polarization assays: 2 kinases, 1 protease, and 1 phosphatase.
Turek-Etienne TC, Small EC, Soh SC, Xin TA, Gaitonde PV, Barrabee EB, Hart RF, Bryant RW
J Biomol Screen (2003) 8:176-176
2. [Patterning enzymes inside microfluidic channels via photoattachment chemistry.](#)
Holden MA, Jung SY, Cremer PS
Anal Chem (2004) 76:1838-1843
3. [Engineered single-chain dimeric streptavidins with an unexpected strong preference for biotin-4-fluorescein.](#)
Aslan FM, Yu Y, Mohr SC, Cantor CR
Proc Natl Acad Sci U S A (2005) 102:8507-8512