

NBD-X, SE (Succinimidyl 6-(N-(7-Nitrobenz-2-Oxa-1,3-Diazol-4-yl)Amino)Hexanoate)

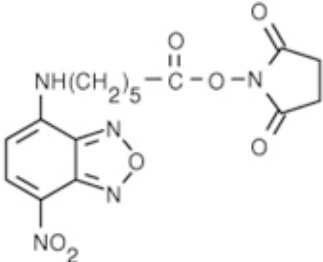
Catalog Number	Packaging Size
C162	25 mg

Storage upon receipt: -20°C, protect from light

Introduction

The amine-reactive **NBD-X, succinimidyl ester** can be used to create environment sensitive bioconjugates. NBD is a functional analog of the dinitrophenyl hapten, and its fluorescence is quenched upon binding to anti-dinitrophenyl (anti-DNP) antibodies.

Specifications

Label:	NBD	
Ex/Em:	466/534 nm	
Detection Method:	Fluorescent	
Solubility:	DMSO, DMF	
Molecular Formula:	C ₁₆ H ₁₇ N ₅ O ₇	
Molecular Weight:	391.34	
CAS Number:	145195-58-0	
Storage Conditions:	-20°C, protect from light	
Shipping Condition:	Room Temperature	

Applications

Fluorescent labeling

References:

1. [Intracellular delivery of phosphoinositides and inositol phosphates using polyamine carriers.](#)
Ozaki S, DeWald DB, Shope JC, Chen J, Prestwich GD
Proc Natl Acad Sci U S A (2000) 97:11286-11291
2. Synthesis and membrane binding properties of a lipopeptide fragment from influenza virus A hemagglutinin.
Eisele F, Kuhlmann J, Waldmann H
Chem Eur J (2002) 8:3362-3362
3. [Solution conformation of a nitrobenzoxadiazole derivative of the polyene antibiotic nystatin: a FRET study.](#)
Silva L, Coutinho A, Fedorov A, Prieto M
J Photochem Photobiol B (2003) 72:17-26