

6-NBDG

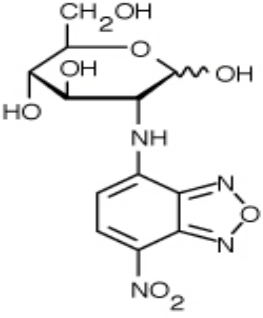
Catalog Number	Packaging Size
C243	5 mg

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

Introduction

6-NBDG is a fluorescent nonhydrolyzable glucose analog that has been used to monitor glucose uptake and transport in live cells. Although sensitive to its environment NBD fluorescence typically displays excitation/emission maxima of ~465/540 nm and can be visualized using optical filters designed for fluorescein.

Specifications

Label:	6-NBDG	
Ex/Em:	465/540 nm	
Detection Method:	Fluorescent	
Solubility:	DMSO, DMF	
Molecular Formula	C ₁₂ H ₁₄ N ₄ O ₈	
Molecular Weight:	342.26	
CAS Number:	108708-22-1	
Storage Conditions:	-20°C, protect from light	
Shipping Condition:	Room Temperature	

Applications

Glucose transport indicator

References:

[Evaluation of glucose transport and its regulation by insulin in human monocytes using flow cytometry.](#)
Dimitriadis G, Maratou E, Boutati E, Psarra K, Papasteriades C, Raptis SA
Cytometry A (2005) 64:27-33

[Glucose Directly Links to Lipid Metabolism through High Affinity Interaction with Peroxisome Proliferator-activated Receptor {alpha}.](#)
Hostetler HA, Huang H, Kier AB, Schroeder F,
J Biol Chem (2008) 283:2246-2254

[Glutamate triggers rapid glucose transport stimulation in astrocytes as evidenced by real-time confocal microscopy.](#)

Loaiza A, Porras OH, Barros LF
J Neurosci (2003) 23:7337-7342