

2-NBD-Glucose

Glucose uptake probe

Fluorescent glucose uptake probe. May be used to monitor glucose uptake in live cells and as a cell viability indicator. Has been used to measure glucose uptake in cultured hippocampal astrocytes using confocal epifluorescence microscopy. NBD fluorescence typically displays excitation/emission maxima of 465/540 nm. Cell permeable.

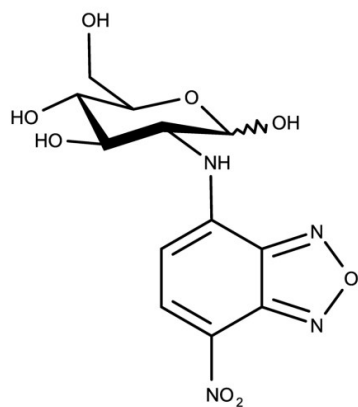
Ordering Information

[Order Online »](#)

ENZ-CHM484-0005	5mg
-----------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Protect from light.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	2-Deoxy-2-[(7-nitro-2,1,3-benzoxadiazol-4-yl)amino]-D-glucose
Appearance	Purple solid.
CAS	186689-07-6
Formula	$C_{12}H_{14}N_4O_8$
Identity	Determined by NMR.
MW	342.27
Purity	≥98% (TLC)
Solubility	Soluble in DMSO (10 mg/ml).

Last modified: February 27, 2025



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-fr@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com