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

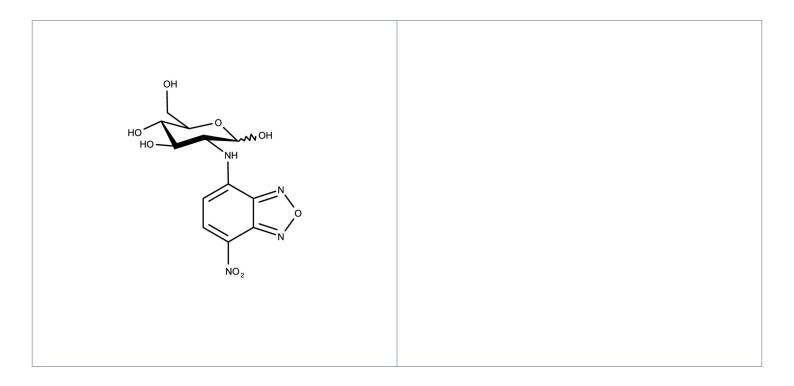
Order Online »

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