

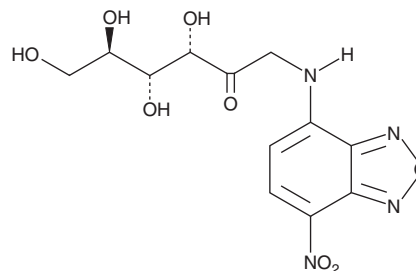
PRODUCT INFORMATION



NBD-Fructose

Item No. 9002314

CAS Registry No.: 940961-04-6
Formal Name: 1-deoxy-1-[(7-nitro-2,1,3-benzoxadiazol-4-yl)amino]-D-fructose
Synonyms: 1-NBDF, 1-(7-nitro-1,2,3-benzadiazole)-Fructose
MF: C₁₂H₁₄N₄O₈
FW: 342.3
Purity: ≥95% (sum of isomers)
UV/Vis.: λ_{max}: 226, 330, 461 nm
Ex./Em. Max: 472/538 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥2 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Description

NBD-Fructose is a fluorescent derivative of fructose that is formed by coupling NBD-chloride with the amine group of amino fructose.¹ It has been used to monitor fructose uptake by the GLUT5 transporter in MCF-7, MDA-MB-435, and MDA-MB-231 human breast cancer cells. NBD-Fructose displays excitation/emission maxima of 472/538 nm, respectively.

Reference

1. Levi, J., Cheng, Z., Gheysens, O., *et al.* Fluorescent fructose derivatives from imaging breast cancer cells. *Bioconjug. Chem.* **18**(3), 628-634 (2007).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 05/30/2018

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM