

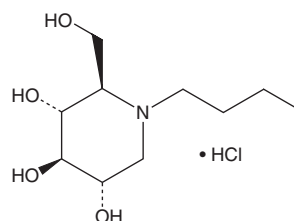
# PRODUCT INFORMATION



## N-Butyldeoxynojirimycin (hydrochloride)

Item No. 21065

**CAS Registry No.:** 210110-90-0  
**Formal Name:** 1-butyl-2R-(hydroxymethyl)-3R,4R,5S-piperidinetriol, monohydrochloride  
**Synonym:** Miglustat, NB-DNJ  
**MF:** C<sub>10</sub>H<sub>21</sub>NO<sub>4</sub> • HCl  
**FW:** 255.7  
**Purity:** ≥98%  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:** ≥4 years



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

### Laboratory Procedures

N-Butyldeoxynojirimycin (NB-DNJ) (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the NB-DNJ (hydrochloride) in the solvent of choice which should be purged with an inert gas, NB-DNJ (hydrochloride) is soluble in the organic solvent DMSO at a concentration of approximately 10 mg/ml.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of NB-DNJ (hydrochloride) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of NB-DNJ (hydrochloride) in PBS, pH 7.2, is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

NB-DNJ is an iminosugar that inhibits UDP-glucose ceramide glucosyltransferase and  $\beta$ -glucosidase 2 ( $IC_{50}$ s = 32 and 81  $\mu$ M, respectively, for the rat recombinant enzymes).<sup>1</sup> It also increases the activity of wild-type acid  $\beta$ -glucosidase, as well as the S364R, N370S, V15M, or M123T mutants, in COS-7 cells expressing the human enzymes when used at a concentration of 10  $\mu$ M.<sup>2</sup> NB-DNJ also inhibits HIV-1 and HIV-2 infection of peripheral blood mononuclear cells (PBMCs;  $IC_{50}$ s = 282 and 211  $\mu$ M, respectively).<sup>3</sup> Formulations containing NB-DNJ have been used in the treatment of Gaucher disease, an inborn error of metabolism characterized as a lysosomal storage disorder resulting from substantial deficiency of  $\beta$ -glucosidase.

### References

1. Lee, J.C., Francis, S., Dutta, D., *et al.* Synthesis and evaluation of eight- and four-membered iminosugar analogues as inhibitors of testicular ceramide-specific glucosyltransferase, testicular  $\beta$ -glucosidase 2, and other glycosidases. *J. Org. Chem.* **77**(7), 3082-3098 (2012).
2. Alfonso, P., Pampín, S., Estrada, J., *et al.* Miglustat (NB-DNJ) works as a chaperone for mutated acid  $\beta$ -glucosidase in cells transfected with several Gaucher disease mutations. *Blood Cells Mol. Dis.* **35**(2), 268-276 (2005).
3. Pollock, S., Dwek, R.A., Burton, D.R., *et al.* N-Butyldeoxynojirimycin is a broadly effective anti-HIV therapy significantly enhanced by targeted liposome delivery. *AIDS* **22**(15), 1961-1969 (2008).

#### 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, 10/07/2022

#### 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](http://WWW.CAYMANCHEM.COM)