

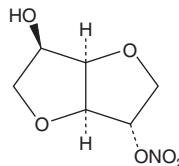
PRODUCT INFORMATION



Isosorbide-2-nitrate

Item No. 9003732

CAS Registry No.: 16106-20-0
Formal Name: 1,4:3,6-dianhydro-D-glucitol, 2-nitrate
Synonyms: IS 2-MN, Isosorbide 2-mononitrate
MF: $C_6H_9NO_6$
FW: 191.1
Purity: $\geq 98\%$
Supplied as: A solid
Storage: -20°C
Stability: ≥ 2 years



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

Laboratory Procedures

Isosorbide-2-nitrate is supplied as a solid. A stock solution may be made by dissolving the isosorbide-2-nitrate in the solvent of choice, which should be purged with an inert gas. Isosorbide-2-nitrate is soluble in methanol and chloroform.

Description

Isosorbide-2-nitrate is a nitric oxide (NO) donor and an active metabolite of isosorbide dinitrate (Item No. 23990).^{1,2} It inhibits potassium chloride- or angiotensin II-induced contraction of rabbit aortic strips and vasopressin-induced coronary spasm in isolated rabbit hearts when used at concentrations of 100 and 1,000 μM .

References

1. Raddino, R., Caretta, G., Bonadei, I., *et al.* Differences between nitrates: Role of isosorbide 2-mononitrate. *J. Cardiovasc. Med. (Hagerstown)* **11**(7), 486-492 (2010).
2. Patel, B.D., Pulakundam, N., Gadekar, A.R., *et al.* Development and validation of a chiral LC-ESI-MS/MS method for simultaneous determination of the enantiomeric metabolites isosorbide 2-mononitrate and isosorbide 5-mononitrate of isosorbide dinitrate in rat and human plasma. *J. Pharm. Biomed. Anal.* **174**, 396-405 (2019).

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, 04/25/2025

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