

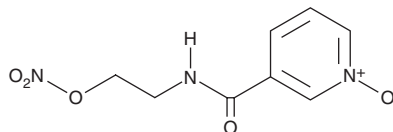
PRODUCT INFORMATION



Nicorandil N-oxide

Item No. 35415

CAS Registry No.: 107833-98-7
Formal Name: N-[2-(nitrooxy)ethyl]-3-pyridinecarboxamide 1-oxide
MF: $C_8H_9N_3O_5$
FW: 227.2
Purity: $\geq 98\%$
UV/Vis.: λ_{max} : 221, 271 nm
Supplied as: A 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

Nicorandil N-oxide is supplied as a solid. A stock solution may be made by dissolving the nicorandil N-oxide in the solvent of choice, which should be purged with an inert gas. Nicorandil N-oxide is slightly soluble in DMSO and methanol.

Description

Nicorandil N-oxide is a metabolite of the sulfonylurea receptor 2B (SUR2B) linked to ATP-sensitive potassium channel $K_{ir}6.2$ (SUR2B/ $K_{ir}6.2$) activator and nitric oxide donor nicorandil (Item No. 18460).¹

Reference

1. Frydman, A.M., Chapelle, P., Diekmann, H., et al. Pharmacokinetics of nicorandil. *Am. J. Cardiol.* **63**(21), 25J-33J (1989).

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, 11/19/2024

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