

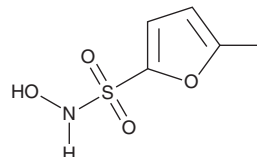
PRODUCT INFORMATION



Cimlanod

Item No. 40660

CAS Registry No.: 1620330-72-4
Formal Name: N-hydroxy-5-methyl-2-furansulfonamide
Synonyms: BMS-986231, CXL-1427
MF: $C_5H_7NO_4S$
FW: 177.2
Purity: $\geq 98\%$
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

Cimlanod is supplied as a solid. A stock solution may be made by dissolving the cimlanod in the solvent of choice, which should be purged with an inert gas. Cimlanod is slightly soluble (0.1-1 mg/ml) in acetonitrile and sparingly soluble (1-10 mg/ml) in DMSO.

Description

Cimlanod is a nitroxyl donor.¹ *In vivo*, cimlanod (7 $\mu\text{g}/\text{kg}/\text{min}$ infusion for four hours) reduces mean arterial pressure (MAP) and systemic vascular resistance and increases cardiac output at four hours post infusion in a dog model of microembolism-induced heart failure.

Reference

1. Hartman, J.C., del Rio, C.L., Reardon, J.E., *et al.* Intravenous infusion of the novel HNO donor BMS-986231 is associated with beneficial inotropic, lusitropic, and vasodilatory properties in 2 canine models of heart failure. *JACC Basic Transl. Sci.* **3**(5), 625-638 (2018).

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, 06/07/2024

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