

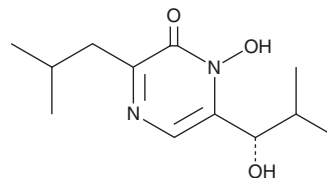
# PRODUCT INFORMATION



## Neohydroxyaspergillic Acid

Item No. 24918

CAS Registry No.: 72598-34-6  
Formal Name: 1-hydroxy-6-[(1S)-1-hydroxy-2-methylpropyl]-3-(2-methylpropyl)-2(1H)-pyrazinone  
Synonym: NSC 613947  
MF:  $C_{12}H_{20}N_2O_3$   
FW: 240.3  
Purity:  $\geq 95\%$   
Supplied as: A solid  
Storage:  $-20^{\circ}\text{C}$   
Stability:  $\geq 4$  years  
Item Origin: *Aspergillus* sp.



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

### Laboratory Procedures

Neohydroxyaspergillic acid is supplied as a solid. A stock solution may be made by dissolving the neohydroxyaspergillic acid in the solvent of choice. Neohydroxyaspergillic acid is soluble in organic solvents such as ethanol, methanol, DMSO, and dimethyl formamide, which should be purged with an inert gas.

### Description

Neohydroxyaspergillic acid is a fungal metabolite produced by *A. sclerotiorum* that has antibiotic and antifungal activities.<sup>1</sup> It inhibits the growth of *P. aeruginosa*, *M. smegmatis*, *S. aureus*, *E. coli*, *K. pneumoniae*, *B. mycoides*, and *B. subtilis* bacteria (MICs = 125-500  $\mu\text{g/ml}$ ). Neohydroxyaspergillic acid also inhibits the growth of *G. convolute*, *S. consortiale*, *P. blakesleeana*, *C. globosum*, and *T. mentagrophytes* (MICs = 175-700  $\mu\text{g/ml}$ ) but not *A. niger*, *P. notatum*, *M. verrucaria*, or *S. cerevisiae* fungi.

### Reference

1. Weiss, U., Strelitz, F., Flon, H., et al. Antibiotic compounds with action against bacterial viruses: Neohydroxyaspergillic acid. *Arch. Biochem. Biophys.* **74**(1), 150-157 (1958).

#### 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, 12/06/2022

#### 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