

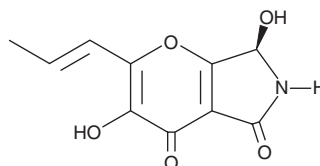
PRODUCT INFORMATION



Pyranonigrin A

Item No. 27535

CAS Registry No.: 773855-65-5
Formal Name: (7R)-6,7-dihydro-3,7-dihydroxy-2-(1E)-1-propen-1-yl-pyrano[2,3-c]pyrrole-4,5-dione
Synonym: 7-hydroxy Pyranonigrin S
MF: C₁₀H₉NO₅
FW: 223.2
Purity: ≥95%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Fungus/*Aspergillus niger*



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

Laboratory Procedures

Pyranonigrin A is supplied as a solid. A stock solution may be made by dissolving the pyranonigrin A in the solvent of choice, which should be purged with an inert gas. Pyranonigrin A is soluble in DMSO and methanol (warmed).

Description

Pyranonigrin A is a fungal metabolite originally isolated from *Aspergillus* that has antioxidant activity.¹ It scavenges 2,2-diphenyl-1-picrylhydrazyl (DPPH; Item No. 14805) free radicals in a cell-free assay (IC₅₀ = 132.9 μM). Pyranonigrin A (10 μM) suppresses TNF-α-induced expression of vascular cell adhesion molecule 1 (VCAM-1) in human umbilical vein endothelial cells (HUVECs).

Reference

1. Miyake, Y., Mochizuki, M., Ito, C., *et al.* Antioxidative pyranonigrins in rice mold starters and their suppressive effect on the expression of blood adhesion molecules. *Biosci. Biotech. Biochem.* **72(6)**, 1580-1585 (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, 08/09/2023

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