

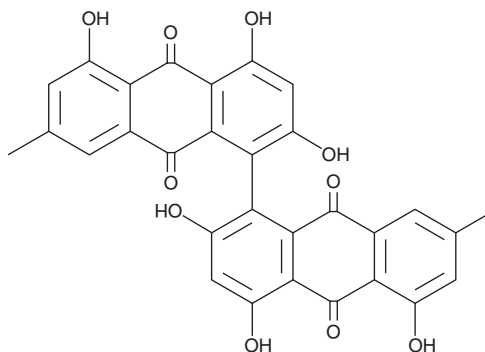
PRODUCT INFORMATION



Skyrin

Item No. 13102

CAS Registry No.: 602-06-2
Formal Name: 2,2',4,4',5,5'-hexahydroxy-7,7'-dimethyl-[1S,1'-bianthracene]-9,9',10,10'-tetrone
Synonyms: Endothianin, Rhodophyscin
MF: C₃₀H₁₈O₁₀
FW: 538.5
Purity: ≥97%
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Fungus/*Penicillium* sp.



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

Laboratory Procedures

Skyrin is supplied as a crystalline solid. A stock solution may be made by dissolving the skyrin in the solvent of choice. Skyrin is soluble in organic solvents such as DMSO, which should be purged with an inert gas.

Description

Skyrin is a fungal metabolite characterized by a bisanthraquinone structure. It interferes with glucagon signaling through adenylate cyclase without binding to the glucagon receptor.¹ At 10 μM, skyrin reduces both glucagon-stimulated cAMP production and glycogenolysis.¹ It does not interfere with epinephrine or glucagon-like peptide 1 effects on these parameters.¹

Reference

1. Parker, J.C., McPherson, R.K., Andrews, K.M., *et al.* Effects of skyrin, a receptor-selective glucagon antagonist, in rat and human hepatocytes. *Diabetes* **49**, 2079-2086 (2012).

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, 08/10/2023

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