

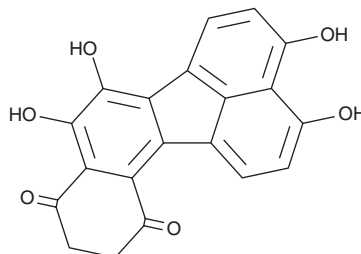
# PRODUCT INFORMATION



## Hortein

Item No. 33250

CAS Registry No.: 346610-88-6  
Formal Name: 2,3-dihydro-5,6,9,10-tetrahydroxy-benzo[j]  
fluoranthene-1,4-dione  
MF:  $C_{20}H_{12}O_6$   
FW: 348.3  
Purity:  $\geq 95\%$   
UV/Vis.:  $\lambda_{\max}$ : 257, 357 nm  
Supplied as: A solid  
Storage:  $-20^{\circ}\text{C}$   
Stability:  $\geq 4$  years  
Item Origin: Fungus/Unidentified sp.



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

### Laboratory Procedures

Hortein is supplied as a solid. A stock solution may be made by dissolving the hortein in the solvent of choice, which should be purged with an inert gas. Hortein is soluble in DMSO.

### Description

Hortein is a polyketide fungal metabolite originally isolated from *H. werneckii*.<sup>1</sup>

### Reference

1. Brauers, G., Ebel, R., Edrada, R., *et al.* Hortein, a new natural product from the fungus *Hortaea werneckii* associated with the sponge *Aplysina aerophoba*. *J. Nat. Prod.* **64**(5), 651-652 (2001).

#### 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, 08/25/2023

#### 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