

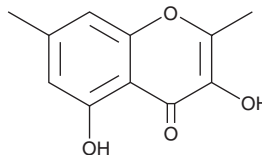
# PRODUCT INFORMATION



## Lachnone A

Item No. 26462

CAS Registry No.: 903892-99-9  
Formal Name: 3,5-dihydroxy-2,7-dimethyl-4H-1-benzopyran-4-one  
MF: C<sub>11</sub>H<sub>10</sub>O<sub>4</sub>  
FW: 206.2  
Purity: ≥95%  
Supplied as: A solid  
Storage: -20°C  
Stability: ≥4 years  
Item Origin: Fungi



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

### Laboratory Procedures

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

Lachnone A is sparingly soluble in aqueous solutions. To enhance aqueous solubility, dilute the organic solvent solution into aqueous buffers or isotonic saline. If performing biological experiments, ensure the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

### Description

Lachnone A is a chromone fungal metabolite originally isolated from *Lachnum*.<sup>1</sup> It is inactive against *P. falciparum*, *C. albicans*, and *M. tuberculosis*. It is not cytotoxic to KB or BC-1 cells.

### Reference

1. Rukachaisirikul, V., Chantaruk, S., Pongcharoen, W., et al. Chromone derivatives from the filamentous fungus *Lachnum* sp. BCC 2424. *J. Nat. Prod.* **69**(6), 980-982 (2006).

#### 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, 12/20/2022

#### 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