

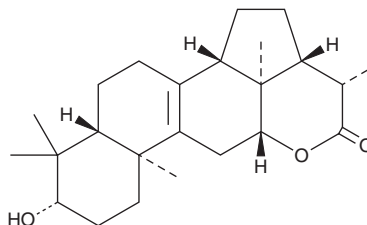
# PRODUCT INFORMATION



## Carbolactone

Item No. 29838

**CAS Registry No.:** 155443-55-3  
**Formal Name:** (5 $\alpha$ )-3 $\beta$ ,12 $\beta$ -dihydroxy-4,4-dimethyl-pregn-8-ene-20S-carboxylic acid,  $\delta$ -lactone  
**MF:** C<sub>24</sub>H<sub>36</sub>O<sub>3</sub>  
**FW:** 372.5  
**Purity:**  $\geq$ 70%  
**Supplied as:** A solid  
**Storage:** -20°C  
**Stability:**  $\geq$ 4 years  
**Item Origin:** Fungus/*Preussia* sp.



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

### Laboratory Procedures

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

### Description

Carbolactone is a fungal metabolite originally isolated from *Sporormiella australis* that has fungicidal and herbicidal activities.<sup>1</sup> It is active against the phytopathogenic fungi *U. violacea* and *E. repens*, but not the bacteria *B. megaterium* or *E. coli*, when used at concentrations of 6.5, 65, and 650  $\mu$ g/ml in an agar diffusion assay. Carbolactone also inhibits germination of the cress *L. sativum*.

### Reference

1. Krohn, K., Ludewig, K., Jones, P.G., *et al.* Biologically active metabolites from fungi, 2<sup>1</sup> an antifungal and herbicidal lanostane lactone from *Sporormiella australis*. *Nat. Prod. Lett.* **1(1)**, 29-32 (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, 08/16/2023

#### 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