

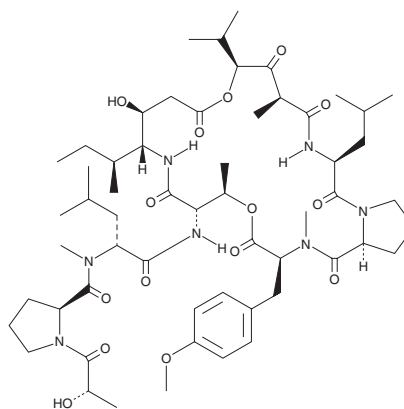
# PRODUCT INFORMATION



## Didemnin B

Item No. 34063

**CAS Registry No.:** 77327-05-0  
**Formal Name:** N-[1-[(2S)-2-hydroxy-1-oxopropyl]-L-prolyl]-didemnin A  
**Synonyms:** NSC 325319, NSC 333841  
**MF:** C<sub>57</sub>H<sub>89</sub>N<sub>7</sub>O<sub>15</sub>  
**FW:** 1,112.4  
**Purity:** ≥98%  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:** ≥1 year  
**Item Origin:** Bacterium/*Tistrella mobilis*



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

### Laboratory Procedures

Didemnin B is supplied as a crystalline solid. A stock solution may be made by dissolving the didemnin B in the solvent of choice, which should be purged with an inert gas. Didemnin B is soluble in the organic solvent acetonitrile at a concentration of approximately 1 mg/ml.

### Description

Didemnin B is a depsipeptide that has been found in sea squirts and has diverse biological activities.<sup>1,2</sup> It reduces viral titers of coxsackievirus A21, equine rhinovirus, parainfluenza virus 3, herpes simplex virus 1 (HSV-1), and HSV-2 in infected Vero cells when used at a concentration of 50 µg/ml.<sup>1</sup> Didemnin B is cytotoxic to murine L1210 lymphocytic leukemia cells (IC<sub>50</sub> = 1.1 ng/ml) and increases survival time in a P388 murine leukemia model when administered at a dose of 0.06 mg/kg. It also reduces splenomegaly in a mouse model of graft versus host reaction (GVHR) when administered at doses ranging from 0.05 to 0.3 mg/kg.<sup>2</sup>

### References

1. Rinehart, K.L., Jr., Gloer, J.B., Hughes, R.G., Jr., *et al.* Didemnins: Antiviral and antitumor depsipeptides from a Caribbean tunicate. *Science* **212**(4497), 933-935 (1981).
2. Montgomery, D.W., and Zukoski, C.L. Didemnin B: A new immunosuppressive cyclic peptide with potent activity in vitro and in vivo. *Transplantation* **40**(1), 49-56 (1985).

#### 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, 08/13/2021

#### 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