

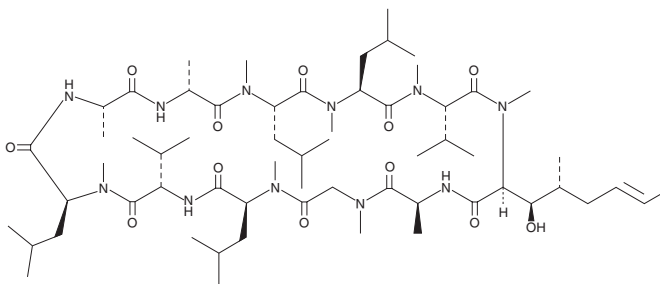
PRODUCT INFORMATION



Cyclosporin B

Item No. 19525

CAS Registry No.: 63775-95-1
Formal Name: 7-L-alanine-cyclosporin A
Synonym: Antibiotic S 7481F2
MF: C₆₁H₁₀₉N₁₁O₁₂
FW: 1,188.6
Purity: ≥98%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

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

Description

Cyclosporin B is a minor cyclopeptide metabolite produced by *T. polysporum* and other deuteromycota with diverse biological activities.¹ It inhibits the growth of T lymphoma cells in a concentration-dependent manner.² Cyclosporin B also inhibits P-glycoprotein (P-gp) activity *in vitro* with an IC₅₀ value of 4.7 μM.³

References

1. Traber, R., Kuhn, M., Loosli, H.R., *et al.* New cyclopeptides from *Trichoderma polysporum* (Link ex Pers.) Rifai: Cyclosporins B, D and E. *Helv. Chim. Acta.* **60(5)**, 1568-1578 (1977).
2. Koponen, M. and Loor, F. Activity of some natural and modified (fluorescent) cyclosporins on T cells. *Ann. Immunol. (Paris)* **134C(2)**, 171-189 (1983).
3. Loor, F., Tiberghien, F., Wenandy, T., *et al.* Cyclosporins: Structure-activity relationships for the inhibition of the human MDR1 P-glycoprotein ABC transporter. *J. Med. Chem.* **45(21)**, 4598-4612 (2002).

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, 08/28/2023

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