

PRODUCT INFORMATION

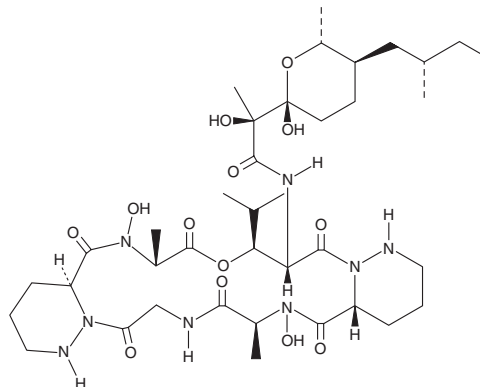


L-156,602

Item No. 23896

CAS Registry No.: 125228-51-5
Formal Name: (3S)-3-hydroxy-N-[(2S)-2-hydroxy-1-oxo-2-[(2R,5R,6R)-tetrahydro-2-hydroxy-6-methyl-5-[(2S)-2-methylbutyl]-2H-pyran-2-yl]propyl]-L-leucyl-(3R)-hexahydro-3-pyridazinocarbonyl-N-hydroxy-L-alanylglycyl-(3S)-hexahydro-3-pyridazinocarbonyl-N-hydroxy-D-alanine, (6→1³)-lactone

Synonym: PD 124966
MF: C₃₈H₆₄N₈O₁₃
FW: 841.0
Purity: ≥95%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Bacterium/*Streptomyces* sp.



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

Laboratory Procedures

L-156,602 is supplied as a solid. A stock solution may be made by dissolving the L-156,602 in the solvent of choice. L-156,602 is soluble in organic solvents such as ethanol, methanol, DMSO, and dimethyl formamide, which should be purged with an inert gas.

Description

L-156,602 is an antagonist of the complement component 5A (C5a) receptor ($IC_{50} = 1.7 \mu\text{g/L}$ in a radioligand binding assay using neutrophil membranes).¹ It is also a cyclic depsipeptide antibiotic that is active against Gram-positive bacteria.² L-156,602 has anti-inflammatory activity in mice, decreasing inflammation induced by muramyl dipeptide (MDP) when administered at a dose of 0.5 mg/kg and by concanavalin A (Item No. 14951) at doses of 0.25 and 0.5 mg/kg.³

References

1. Durette, P.L., Kopka, I.E., Lanza, T.J., *et al.* The inhibition of C5a receptor binding by analogs of L-156,602, a cyclic hexadepsipeptide antibiotic. *Bioor. Med. Chem. Lett.* **2(9)**, 1033-1038 (1992).
2. Hensens, O.D., Borris, R.P., Koupal, L.R., *et al.* L-156,602, a C5a antagonist with a novel cyclic hexadepsipeptide structure from *Streptomyces* sp. MA6348. Fermentation, isolation and structure determination. *J. Antibiot. (Tokyo)* **44(2)**, 249-254 (1991).
3. Tsuji, R.F., Magae, J., Nagai, K., *et al.* Effects of L-156,602, a C5a receptor antagonist, on mouse experimental models of inflammation. *Biosci. Biotech. Biochem.* **56(12)**, 2034-2036 (1992).

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, 10/31/2022

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