

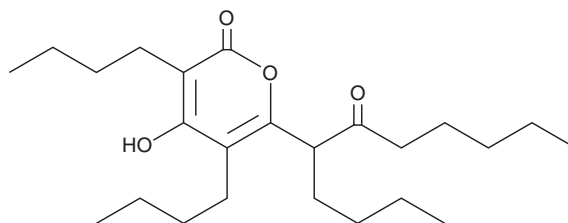
PRODUCT INFORMATION



Elasnin

Item No. 21481

CAS Registry No.: 68112-21-0
Formal Name: 3,5-dibutyl-6-(1-butyl-2-oxoheptyl)-4-hydroxy-2H-pyran-2-one
MF: $C_{24}H_{40}O_4$
FW: 392.6
Purity: $\geq 95\%$
Supplied as: A liquid
Storage: -20°C
Stability: ≥ 2 years
Special Conditions: Light sensitive
Item Origin: Bacterium/*Streptomyces* sp.



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

Laboratory Procedures

Elasnin is supplied as a liquid. A stock solution may be made by dissolving the elasnin in an organic solvent purged with an inert gas. Elasnin is soluble in organic solvents such as methanol and DMSO. The solubility of elasnin in these solvents is approximately 5 mg/ml. Elasnin is also soluble in methylene chloride.

Description

Elasnin is a product isolated from certain strains of *Streptomyces* that acts as a reversible inhibitor of elastase with IC_{50} values of 1.3 and 30 $\mu\text{g/ml}$ for human granulocyte and pancreatic enzymes, respectively.¹ It inhibits human leukocyte elastase, porcine pancreatic elastase, and chymotrypsin with K_i values of 93, 69, and 224 μM , respectively.²

References

1. Ohno, H., Saheki, T., Awaya, J., *et al.* Isolation and characterization of elasnin, a new human granulocyte elastase inhibitor produced by a strain of *Streptomyces*. *J. Antibiot. (Tokyo)* **31(11)**, 1116-1123 (1978).
2. Spencer, R.W., Copp, L.J., and Pfister, J.R. Inhibition of human leukocyte elastase, porcine pancreatic elastase, and chymotrypsin by elasnin and other 4-hydroxy-2-pyrones. *J. Med. Chem.* **28(12)**, 1828-1832 (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, 05/03/2022

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