

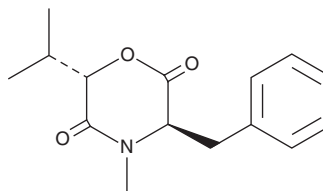
PRODUCT INFORMATION



(3R,6S)-Bassiatin

Item No. 43302

CAS Registry No.: 172721-96-9
Formal Name: 4-methyl-6S-(1-methylethyl)-3R-(phenylmethyl)-2,5-morpholinedione
MF: $C_{15}H_{19}NO_3$
FW: 261.3
Purity: $\geq 98\%$
Supplied as: A solid
Storage: $-20^{\circ}C$
Stability: ≥ 4 years



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

Laboratory Procedures

(3R,6S)-Bassiatin is supplied as a solid. A stock solution may be made by dissolving the (3R,6S)-bassiatin in the solvent of choice, which should be purged with an inert gas. (3R,6S)-Bassiatin is soluble (≥ 10 mg/ml) in DMSO and sparingly soluble (1-10 mg/ml) in ethanol.

Description

(3R,6S)-Bassiatin is the enantiomer of the fungal metabolite bassiatin.^{1,2}

References

1. Kagamizono, T., Nishino, E., Matsumoto, K., *et al.* Bassiatin, a new platelet aggregation inhibitor produced by *Beauveria bassiana* K-717. *J. Antibiot. (Tokyo)* **48(12)**, 1407-1412 (1995).
2. Ola, A.R.B., Aly, A.H., Lin, W., *et al.* Structural revision and absolute configuration of lateritin. *Tetrahedron Lett.* **55(45)**, 6184-6187 (2014).

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, 05/16/2025

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