

PRODUCT INFORMATION



Auristatin PE

Item No. 45983

CAS Registry No.: 149606-27-9

Formal Name: N,N-dimethyl-L-valyl-N-[(1S,2R)-2-methoxy-4-[(2S)-2-[(1R,2R)-1-methoxy-2-methyl-3-oxo-3-[(2-phenylethyl)amino] propyl]-1-pyrrolidinyl]-1-[(1S)-1-methylpropyl]-4-oxobutyl]-N-methyl-L-valinamide

Synonyms: NSC 654663, Soblidotin, TZT-1027, YHI-501

MF: $C_{39}H_{67}N_5O_6$

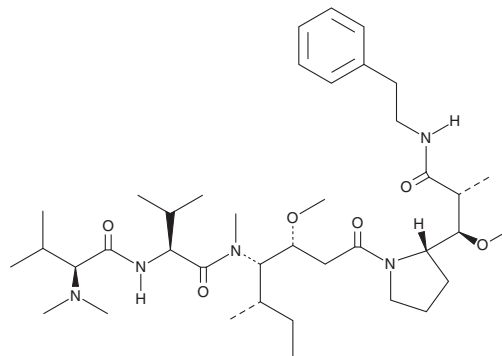
FW: 702.0

Purity: $\geq 98\%$

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

Auristatin PE is supplied as a solid. A stock solution may be made by dissolving the auristatin PE in the solvent of choice, which should be purged with an inert gas. Auristatin PE is slightly soluble (0.1-1 mg/ml) in acetonitrile and DMSO.

Description

Auristatin PE is a tubulin polymerization inhibitor, an anticancer agent, and a derivative of dolastatin 10.¹ It inhibits tubulin polymerization ($IC_{50} < 2.2 \mu\text{M}$), completely inhibits the proliferation of WSU-WM Waldenström's macroglobulinemia cells at 100 pg/ml, and induces cell cycle arrest at the G₂/M phase and apoptosis in the same cells at 10 pg/ml.^{1,2} Auristatin PE induces tumor regression in MX-1 breast cancer, LX-1 lung cancer, and WSU-WM mouse xenograft models.

References

1. Kobayashi, M., Natsume, T., Tamaoki, S., *et al.* Antitumor activity of TZT-1027, a novel dolastatin 10 derivative. *Jpn. J. Cancer Res.* **88**(3), 316-327 (1997).
2. Mohammad, R.M., Limvarapuss, C., Wall, N.R., *et al.* A new tubulin polymerization inhibitor, auristatin PE, induces tumor regression in a human Waldenström's macroglobulinemia xenograft model. *Int. J. Oncol.* **15**, 367-372 (1999).

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, 07/28/2026

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