

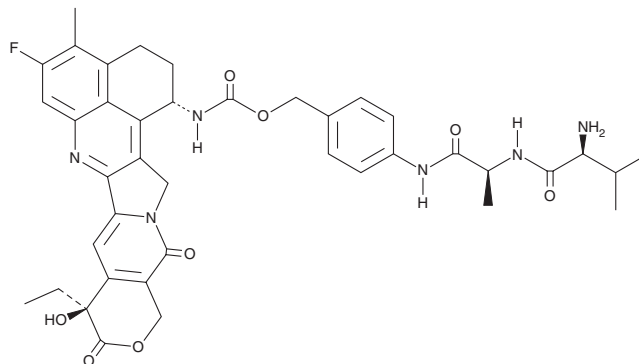
PRODUCT INFORMATION



Val-Ala-PABC-Exatecan

Item No. 44443

CAS Registry No.: 2845164-91-0
Formal Name: L-valyl-N-[4-[[[(1S,9S)-9-ethyl-5-fluoro-2,3,9,10,13,15-hexahydro-9-hydroxy-4-methyl-10,13-dioxo-1H,12H-benzo[de]pyrano[3',4':6,7]indolizino[1,2-b]quinolin-1-yl]amino]carbonyl]oxy]methyl]phenyl]-L-alaninamide
MF: C₄₀H₄₃FN₆O₈
FW: 754.8
Purity: ≥95%
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

Val-Ala-PABC-Exatecan is supplied as a solid. A stock solution may be made by dissolving the Val-Ala-PABC-exatecan in the solvent of choice, which should be purged with an inert gas. Val-Ala-PABC-Exatecan is soluble (≥10 mg/ml) in DMSO.

Description

Val-Ala-PABC-Exatecan is a drug-linker molecule containing the DNA topoisomerase I inhibitor exatecan (Item No. 35452) and a Val-Ala-*p*-aminobenzyl carbamate (Val-Ala-PABC) linker. It has been used in the synthesis of Mal-PEG₈-Val-Ala-PAB-exatecan (Item No. 44595), which has been incorporated into antibody-drug conjugates (ADCs) with anticancer activity in mice.¹

Reference

1. Li, X., Bai, L., Li, X., *et al.* NH₂-trimethyl lock (NH₂-TML): A stable and effective adapter for exatecan-based antibody-drug conjugates. *Bioorg. Chem.* **179**, 110038 (2026).

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, 07/30/2026

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