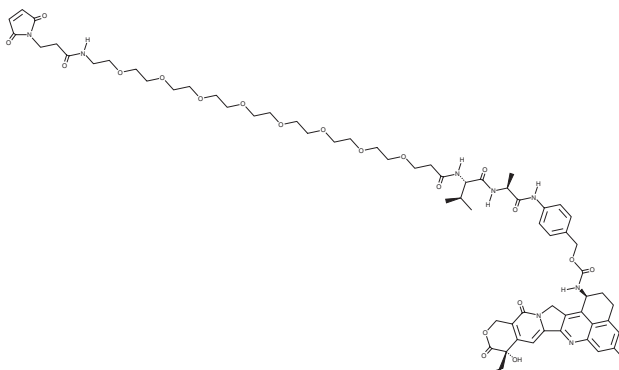


PRODUCT INFORMATION



Mal-PEG₈-Val-Ala-PAB-Exatecan Item No. 44595

CAS Registry No.: 2679821-39-5
Formal Name: 4-((2S,5S)-37-(2,5-dioxo-2,5-dihydro-1H-pyrrol-1-yl)-5-isopropyl-2-methyl-4,7,35-trioxo-10,13,16,19,22,25,28,31-octaoxa-3,6,34-triazaheptatriacontanamido)benzyl ((1S,9S)-9-ethyl-5-fluoro-9-hydroxy-4-methyl-10,13-dioxo-2,3,9,10,13,15-hexahydro-1H,12H-benzo[de]pyrano[3',4':6,7]indolizino[1,2-b]quinolin-1-yl)carbamate
MF: C₆₆H₈₅FN₈O₂₀
FW: 1,329.4
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

Mal-PEG₈-Val-Ala-PAB-Exatecan is supplied as a solid. A stock solution may be made by dissolving the Mal-PEG₈-Val-Ala-PAB-exatecan in the solvent of choice, which should be purged with an inert gas. Mal-PEG₈-Val-Ala-PAB-Exatecan is slightly soluble (0.1-1 mg/ml) in DMSO and methanol.

Description

Mal-PEG₈-Val-Ala-PAB-Exatecan is a reactive drug-linker molecule containing the DNA topoisomerase I inhibitor exatecan (Item No. 35452) and a maleimide-terminal linker.

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, 01/20/2026

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