

PRODUCT INFORMATION



Deruxtecan

Item No. 43189

CAS Registry No.: 1599440-13-7
Formal Name: N-[6-(2,5-dihydro-2,5-dioxo-1H-pyrrol-1-yl)-1-oxohexyl]glycylglycyl-L-phenylalanyl-N-[[2-[[[(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]-2-oxoethoxy)methyl]-glycinamide

Synonyms: Maleimidocaproyl-Gly-Gly-Phe-Gly-Dxd, Mc-GGFG-Dxd

MF: C₅₂H₅₆FN₉O₁₃

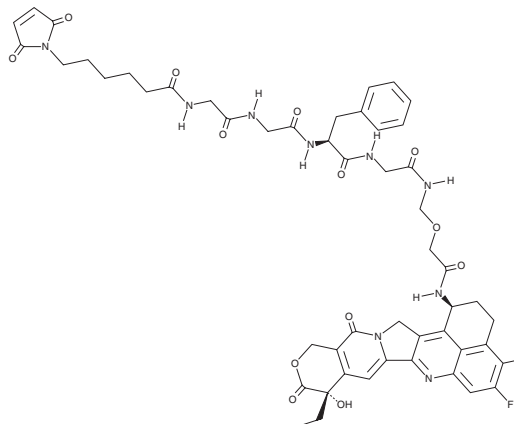
FW: 1,034.1

Purity: ≥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

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

Description

Deruxtecan is a reactive drug-linker molecule containing the exatecan derivative Dxd (Item No. 41699) and a maleimidocaproyl (Mc) linker.¹ It has been linked to the anti-HER2 antibody trastuzumab as an antibody-drug conjugate (ADC) to target Dxd to tumor cells. Deruxtecan linked to trastuzumab binds to HER2 in a cell-free assay, is cytotoxic to breast cancer cells *in vitro*, and induces tumor regression in a BT474 breast cancer mouse xenograft model. Formulations containing deruxtecan conjugated to tumor-targeting antibodies have been used in the treatment of cancer.

Reference

1. Conilh, L., Fournet, G., Fourmaux, E., *et al.* Exatecan antibody drug conjugates based on a hydrophilic polysarcosine drug-linker platform. *Pharmaceuticals (Basel)* **14**(3), 247 (2021).

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, 05/22/2025

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