

PRODUCT INFORMATION



CL2-SN-38 (dichloroacetic acid salt)

Item No. 28215

Formal Name: N-[32-[4-[[[4-[(2,5-dihydro-2,5-dioxo-1H-pyrrol-1-yl)methyl]cyclohexyl]carbonyl]amino]methyl]-1H-1,2,3-triazol-1-yl]-1,5-dioxo-3,9,12,15,18,21,24,27,30-nonaoxa-6-azadotriacont-1-yl]-L-phenylalanyl-N-[4-[[[[(4S)-4,11-diethyl-3,4,12,14-tetrahydro-9-hydroxy-3,14-dioxo-1H-pyrano[3',4':6,7]indolizino[1,2-b]quinolin-4-yl]oxy]carbonyl]oxy]methyl]phenyl]-L-lysineamide, 2,2-dichloro-acetic acid

MF: C₈₂H₁₀₆N₁₂O₂₃ • XC₂H₂Cl₂O₂

FW: 1,627.8

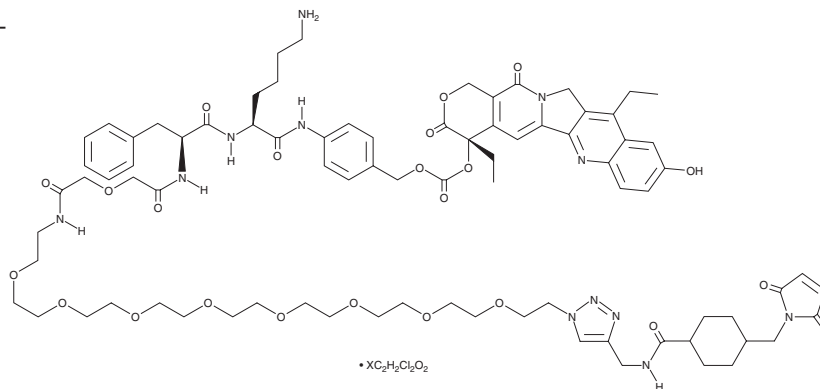
Purity: ≥98%

UV/Vis.: λ_{max}: 223, 268, 329, 367, 382 nm

Supplied as: A crystalline 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

CL2-SN-38 (dichloroacetic acid salt) is supplied as a crystalline solid. A stock solution may be made by dissolving the CL2-SN-38 (dichloroacetic acid salt) in the solvent of choice, which should be purged with an inert gas. CL2-SN-38 (dichloroacetic acid salt) is soluble in organic solvents such as ethanol and DMSO. The solubility of CL2-SN-38 (dichloroacetic acid salt) in these solvents is approximately 1 mg/ml.

Description

CL2-SN-38 is a cleavable linker-based antibody-drug conjugate (ADC) containing the topoisomerase I inhibitor SN-38 (Item No. 15632) and a maleimidocaproyl linker.¹ CL2-SN-38 has been linked to tumor-selective human monoclonal antibodies, such as anti-CEACAM5 (hMN-14) to target SN-38 to tumor sites and improve its solubility. CL2-SN-38 linked to hMN-14 binds (K_d = 1.4 nM) and is cytotoxic to LoVo human colon adenocarcinoma cells (IC₅₀ = 5.3 nM). CL2-SN-38 linked to hMN-14 increases survival in a colon carcinoma mouse model of lung metastasis.

Reference

1. Moon, S., Govindan, S.V., Cardillo, T.M., *et al.* Antibody conjugates of 7-ethyl-10-hydroxycamptothecin (SN-38) for targeted cancer chemotherapy. *J Med Chem.* **51**(21), 6916–6926 (2008).

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, 12/08/2022

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