

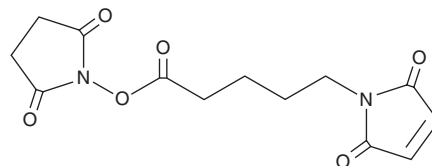
PRODUCT INFORMATION



5-Maleimido-pentanoic NHS ester

Item No. 41900

CAS Registry No.: 103750-03-4
Formal Name: 2,5-dihydro-2,5-dioxo-1H-pyrrole-1-pentanoic acid, 2,5-dioxo-1-pyrrolidinyl ester
Synonyms: 5-Maleimido-pentanoic N-hydroxysuccinimide ester, 5-Maleimido Valeric Acid-NHS, 5-Maleimido Valeric Acid-N-hydroxysuccinimide
MF: C₁₃H₁₄N₂O₆
FW: 294.3
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

5-Maleimido-pentanoic NHS ester is supplied as a solid. A stock solution may be made by dissolving the 5-maleimido-pentanoic NHS ester in the solvent of choice, which should be purged with an inert gas. 5-Maleimido-pentanoic NHS ester is soluble (≥10 mg/ml) in DMSO.

Description

5-Maleimido-pentanoic NHS ester is a linker molecule.¹ It has been used to conjugate SN-38 (Item No. 15632) to phlorizin (Item No. 11576) or phloretin (Item No. 14452) to increase the selectivity of SN-38 for cancer cells.

Reference

1. Chiu, P.F., Chang, C.K., Huang, P.S., *et al.* Design, synthesis, and evaluation of glucose transporter inhibitor-SN38 conjugates for targeting colorectal cancer. *J. Med. Chem.* **66**(14), 9681-9696 (2023).

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, 10/03/2024

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