

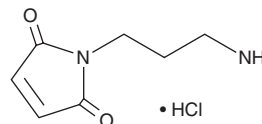
PRODUCT INFORMATION



N-(2-Aminopropyl)maleimide (hydrochloride)

Item No. 45145

CAS Registry No.: 1257852-74-6
Formal Name: 1-(3-aminopropyl)-1H-pyrrole-2,5-dione, monohydrochloride
Synonym: Mal-NH₂
MF: C₇H₁₀N₂O₂ • HCl
FW: 190.6
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

N-(2-Aminopropyl)maleimide (Mal-NH₂) (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the mal-NH₂ (hydrochloride) in the solvent of choice, which should be purged with an inert gas. Mal-NH₂ (hydrochloride) is soluble (≥10 mg/ml) in DMSO.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of mal-NH₂ (hydrochloride) (hydrochloride) can be prepared by directly dissolving the solid in aqueous buffers. Mal-NH₂ (hydrochloride) (hydrochloride) is soluble (≥10 mg/ml) in PBS (pH 7.2). We do not recommend storing the aqueous solution for more than one day.

Description

Mal-NH₂ is a heterobifunctional cross-linking reagent that contains both a thiol-reactive maleimide group and an amine.¹ Mal-NH₂ has been used in the synthesis of polyprodrug nanoparticles as anticancer agents.

Reference

1. Guo, C., Cheng, X., Yang, Y., *et al.* Aptamer-modified GSH-degradable honokiol polyprodrug nanoparticles for ovarian cancer-specific targeting therapy. *Bioorg. Med. Chem. Lett.* **123**, 130215 (2025).

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, 02/11/2026

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