

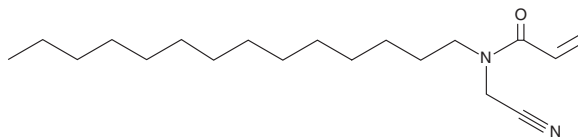
PRODUCT INFORMATION



Cyano-myrcrylamide

Item No. 35155

CAS Registry No.: 2801702-34-9
Formal Name: N-(cyanomethyl)-N-tetradecyl-2-propenamamide
Synonym: CMA
MF: C₁₉H₃₄N₂O
FW: 306.5
Purity: ≥98%
UV/Vis.: λ_{max}: 234 nm
Supplied as: A solid
Storage: -20°C
Stability: ≥2 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

Cyano-myrcrylamide (CMA) is supplied as a solid. A stock solution may be made by dissolving the CMA in the solvent of choice, which should be purged with an inert gas. CMA is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of CMA in these solvents is approximately 30 mg/ml.

Description

CMA is an inhibitor of zinc finger DHHC domain-containing (zDHHC) palmitoyltransferase 20 (zDHHC20; IC₅₀ = 1.35 μM in a cell-free assay).¹ It selectively inhibits zDHHC20 over acyl-protein thioesterase 1 (APT-1) and APT-2 at 50 μM. CMA (5-20 μM) inhibits S-acylation of *Legionella* E3 ligase GobX, MyD88, and Ras, which are substrates of zDHHC20, zDHHC9, and zDHHC6, respectively, in HEK293T cells expressing recombinant *Legionella* GobX, recombinant human MyD88, or endogenous Ras. Unlike the zDHHC inhibitor 2-bromopalmitate (2BP), CMA is not toxic to HEK293T cells at 40 μM.

Reference

1. Azizi, S.-A., Lan, T., Delalande, C., *et al.* Development of an acrylamide-based inhibitor of protein S-acylation. *ACS Chem. Biol.* **16**(8), 1546-1556 (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, 03/30/2023

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