

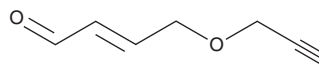
PRODUCT INFORMATION



yne-ACR

Item No. 43702

CAS Registry No.: 220899-12-7
Formal Name: (2E)-4-(2-propyn-1-yloxy)-2-butenal
Synonym: yne-Acrolein
MF: $C_7H_8O_2$
FW: 124.1
Purity: $\geq 98\%$
Supplied as: A 10 mg/ml solution in methyl acetate
Storage: -20°C
Stability: ≥ 2 years



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

Description

yne-ACR is a probe for simulating lipid-derived electrophiles (LDEs) that contains a clickable tag and an electrophilic acrolein warhead.¹ It has been used to modify interfacial and non-adsorbed proteins in a model of interface-dominated systems (IDSs) to study protein behavior during and after LDE-mediated modification.

Reference

1. Gong, W., Jiang, K., Yang, T.-Q., *et al.* Lipid-derived electrophiles modify proteins and alter their interfacial behavior: The distinct mediating role of the interface. *ACS Nano* **19**(16), 15596-15616 (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, 09/22/2025

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