

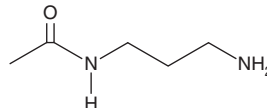
PRODUCT INFORMATION



N-(3-Aminopropyl)acetamide

Item No. 9002787

CAS Registry No.: 4078-13-1
Formal Name: N-(3-aminopropyl)-acetamide
MF: $C_5H_{12}N_2O$
FW: 116.2
Purity: $\geq 95\%$
Supplied as: A solution in ethanol
Storage: $-20^\circ C$
Stability: ≥ 2 years



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

Laboratory Procedures

N-(3-Aminopropyl)acetamide is supplied as a solution in ethanol. To change the solvent, simply evaporate the ethanol under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as ethanol, DMSO, and dimethyl formamide purged with an inert gas can be used. The solubility of N-(3-Aminopropyl)acetamide in these solvents is approximately 30 mg/ml.

Description

N-(3-Aminopropyl)acetamide is a monoacetylated polyamine that can be used as a building block in the synthesis of heterocyclic compounds. Polyamines are promising biochemical markers of cancer and many other pathophysiological conditions, thus their concentrations in biological fluids have been analyzed.¹

Reference

1. Häkkinen, M. R., Roine, A., Auriola, S., *et al.* Analysis of free, mono- and diacetylated polyamines from human urine by LC-MS/MS. *J. Chromatogr. B. Analyt. Technol. Biomed. Life Sci.* (941), 81-89 (2015).

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, 11/11/2022

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