

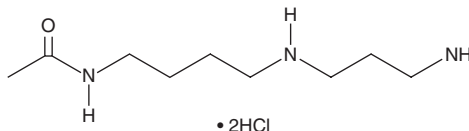
PRODUCT INFORMATION



N⁸- Acetylspermidine (hydrochloride)

Item No. 37434

CAS Registry No.: 34450-15-2
Formal Name: N-[4-[(3-aminopropyl)amino]butyl]-acetamide, dihydrochloride
MF: C₉H₂₁N₃O • 2HCl
FW: 260.2
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

N⁸-Acetylspermidine (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the N⁸-acetylspermidine (hydrochloride) in the solvent of choice, which should be purged with an inert gas. N⁸-Acetylspermidine (hydrochloride) is slightly soluble in methanol and DMSO.

Description

N⁸-Acetylspermidine is an endogenous polyamine.¹⁻³ Urinary levels of N⁸-acetylspermidine are increased in patients with chronic pancreatitis.¹ Blood levels of N⁸-acetylspermidine are increased in patients with ischemic cardiomyopathy, as well as in patients with Parkinson's disease.^{2,3}

References

1. Nissinen, S.I., Venalainen, M., Kumpulainen, P., *et al.* Discrimination between pancreatic cancer, pancreatitis and healthy controls using urinary polyamine panel. *Cancer Control* **29**, 1-8 (2021).
2. Saiki, S., Sasazawa, Y., Fujimaki, M., *et al.* A metabolic profile of polyamines in parkinson disease: A promising biomarker. *Ann. Neurol.* **86**(2), 251-263 (2019).
3. Nayak, A., Liu, C., Mehta, A., *et al.* N8-Acetylspermidine: A polyamine biomarker in ischemic cardiomyopathy with reduced ejection fraction. *J. Am. Heart Assoc.* **9**(11), e016055 (2020).

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, 10/20/2022

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