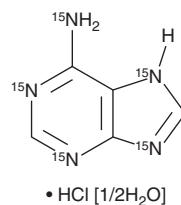


PRODUCT INFORMATION



Adenine-¹⁵N₅ (hydrochloride hydrate) Item No. 41181

Formal Name: 7H-purin-¹⁵N₄-6-amine, monohydrochloride, hemihydrate
MF: C₅H₅[¹⁵N]₅ • HCl [1/2H₂O]
FW: 185.6
Purity: ≥98%
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

Adenine-¹⁵N₅ is intended for use as an internal standard for the quantification of adenine (Item No. 18148) by GC- or LC-MS. Adenine-¹⁵N₅ (hydrochloride hydrate) is supplied as a solid. A stock solution may be made by dissolving the adenine-¹⁵N₅ (hydrochloride hydrate) in the solvent of choice, which should be purged with an inert gas. Adenine-¹⁵N₅ (hydrochloride hydrate) is slightly soluble (0.1-1 mg/ml) in acetonitrile.

Description

Adenine is a purine base that acts as a precursor, substrate, or cofactor in diverse biochemical pathways.¹ It is a fundamental component of DNA, RNA, and ATP, as well as numerous other natural biochemicals.

Reference

1. Nuki, G. Human purine metabolism: Some recent advances and relationships with immunodeficiency. *Ann. Rheum. Dis.* **42 (Suppl 1(Suppl 1))**, 8-11 (1983).

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, 06/25/2024

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