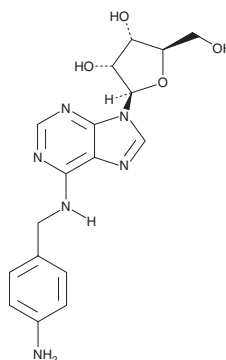


PRODUCT INFORMATION

N-[(4-Aminophenyl)methyl]adenosine

Item No. 28417

CAS Registry No.: 95523-13-0
Synonyms: N-[(4-Aminobenzyl)methyl]adenosine,
 N⁶-4-Aminobenzyladenosine,
 N⁶-p-Aminobenzyladenosine
MF: C₁₇H₂₀N₆O₄
FW: 372.4
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

N-[(4-Aminophenyl)methyl]adenosine is supplied as a solid. A stock solution may be made by dissolving the N-[(4-aminophenyl)methyl]adenosine in the solvent of choice, which should be purged with an inert gas. N-[(4-Aminophenyl)methyl]adenosine is sparingly soluble (1-10 mg/ml) in DMSO.

Description

N-[(4-Aminophenyl)methyl]adenosine is a derivative of adenosine (Item No. 21232) and an intermediate in the synthesis of the radioligand and adenosine receptor agonist [¹²⁵I]-N⁶-4-aminobenzyladenosine.¹

Reference

1. Linden, J., Patel, A., and Sadek, S. [¹²⁵I]Aminobenzyladenosine, a new radioligand with improved specific binding to adenosine receptors in heart. *Circ. Res.* **56(2)**, 279-284 (1985).

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, 10/11/2024

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