

PRODUCT INFORMATION

Adenosine-5'-O-triphosphoro-γ-1-(5-sulfonic acid)naphthylamidate (sodium salt)

Item No. 38537

Formal Name: adenosine 5'-(trihydrogen diphosphate)
P'→P-anhydride with 5-(phosphonoamino)-

Synonym:

ATP-γ-Aminonaphthalene-5-Sulfonic Acid

MF: C₂₀H₂₀N₆O₁₅P₃S • 3Na

FW: 778.4

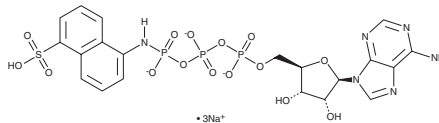
Purity: ≥95%

Ex./Em. Max: 320/460 nm

Supplied as: A solution in water

Storage: -80°C

Stability: ≥4 years



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

Description

Adenosine-5'-O-triphosphoro-γ-1-(5-sulfonic acid)naphthylamidate is a fluorescent derivative of adenosine 5'-triphosphate (Item No. 14498).¹ It displays an emission maximum of 460 nm upon excitation at 320 nm.¹

Reference

1. Yarbrough, L.R. Synthesis and properties of a new fluorescent analog of ATP: Adenosine-5'-triphosphoro-γ-1-(5-sulfonic acid) naphthylamidate. *Biochem. Biophys. Res. Commun.* **81**(1), 35-41 (1978).

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, 04/21/2026

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