

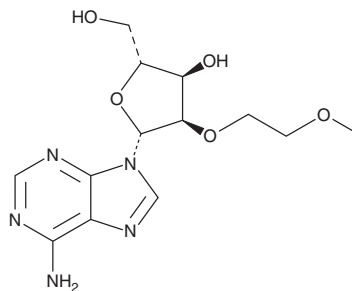
PRODUCT INFORMATION



2'-O-(2-Methoxyethyl)adenosine

Item No. 40443

CAS Registry No.: 168427-74-5
Synonym: 2'-MOE-Adenosine
MF: C₁₃H₁₉N₅O₅
FW: 325.3
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

2'-O-(2-Methoxyethyl)adenosine is supplied as a solid. A stock solution may be made by dissolving the 2'-O-(2-methoxyethyl)adenosine in the solvent of choice, which should be purged with an inert gas. 2'-O-(2-Methoxyethyl)adenosine is soluble in DMSO. 2'-O-(2-Methoxyethyl)adenosine is slightly soluble in acetonitrile.

2'-O-(2-Methoxyethyl)adenosine is slightly soluble in aqueous solutions. To enhance aqueous solubility, dilute the organic solvent solution into aqueous buffers or isotonic saline. If performing biological experiments, ensure the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

Description

2'-O-(2-Methoxyethyl)adenosine is a 2'-O-methoxyethyl derivative of the purine nucleoside adenosine (Item No. 21232).¹ 2'-O-(2-Methoxyethyl)adenosine has been used in VEGF-binding aptamers.

Reference

1. Freund, N., Taylor, A.I., Arangundy-Franklin, S., *et al.* A two-residue nascent-strand steric gate controls synthesis of 2'-O-methyl- and 2'-O-(2-methoxyethyl)-RNA. *Nat. Chem.* **15**(1), 91-100 (2023).

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, 03/08/2024

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