

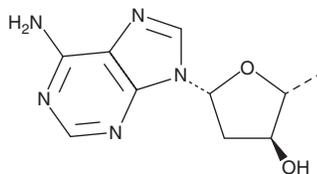
# PRODUCT INFORMATION



## 2',5'-dideoxy Adenosine

Item No. 20358

**CAS Registry No.:** 6698-26-6  
**Formal Name:** 2',5'-dideoxy-adenosine  
**Synonyms:** 2',5'-dd-Ado, NSC 95943  
**MF:** C<sub>10</sub>H<sub>13</sub>N<sub>5</sub>O<sub>2</sub>  
**FW:** 235.2  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 210, 259 nm  
**Supplied as:** A crystalline 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',5'-dideoxy Adenosine is supplied as a crystalline solid. A stock solution may be made by dissolving the 2',5'-dideoxy adenosine in the solvent of choice, which should be purged with an inert gas. 2',5'-dideoxy Adenosine is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of 2',5'-dideoxy adenosine in these solvents is approximately 20 and 5 mg/ml, respectively.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of 2',5'-dideoxy adenosine can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of 2',5'-dideoxy adenosine in PBS, pH 7.2, is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

2',5'-dideoxy Adenosine is a nucleoside analog that is one of the first identified cell-permeable, P-site inhibitors of adenylate cyclase. It inhibits forskolin-induced activation of a cAMP-dependent reporter gene in HEK293 cells with an IC<sub>50</sub> value of 33 μM.<sup>1</sup>

### Reference

1. Emery, A.C., Eiden, M.V., and Eiden, L.E. A new site and mechanism of action for the widely used adenylate cyclase inhibitor SQ22,536. *Mol.Pharmacol.* **83(1)**, 95-105 (2013).

#### 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, 12/01/2022

#### 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