

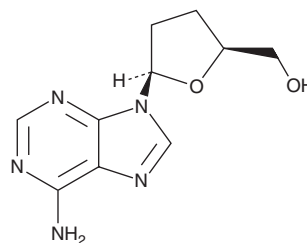
PRODUCT INFORMATION



2',3'-Dideoxyadenosine

Item No. 36662

CAS Registry No.: 4097-22-7
Synonyms: ddA, ddAdo, NSC 98700
MF: $C_{10}H_{13}N_5O_2$
FW: 235.2
Purity: $\geq 98\%$
UV/Vis.: λ_{max} : 262 nm
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',3'-Dideoxyadenosine is supplied as a solid. A stock solution may be made by dissolving the 2',3'-dideoxyadenosine in the solvent of choice, which should be purged with an inert gas. 2',3'-Dideoxyadenosine is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of 2',3'-dideoxyadenosine in these solvents is approximately 2 and 1 mg/ml, respectively. It is slightly soluble in ethanol.

Description

2',3'-Dideoxyadenosine is prodrug form of the nucleoside reverse transcriptase inhibitor (NRTI) 2',3'-dideoxyadenosine 5'-triphosphate (ddATP; Item No. 17460).¹ It inhibits HIV replication in HIV-infected MT-4 cells when used at concentrations of 5 and 125 μM and protects ATH8 cells against the cytopathic effect of HIV ($\text{EC}_{50} = 2.7 \mu\text{M}$).²

References

- Schlienger, N., Lefebvre, I., Aubertin, A.-M., *et al.* Mononucleoside phosphorodithiolates as mononucleotide prodrugs. *Eur. J. Med. Chem.* **227**, 113914 (2022).
- Herdewijn, P., Blazarini, J., Clercq, E.D., *et al.* 3'-substituted 2',3'-dideoxynucleoside analogues as potential anti-HIV (HTLV-III/LAV) agents. *J. Med. Chem.* **30(8)**, 1270-1278 (1987).

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, 12/20/2022

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