

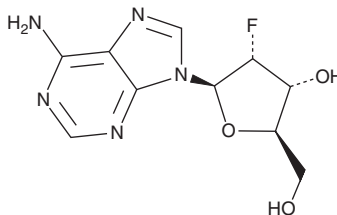
PRODUCT INFORMATION



2'-Fluoro-2'-deoxyadenosine

Item No. 39923

CAS Registry No.: 64183-27-3
Formal Name: 2'-deoxy-2'-fluoro-adenosine
Synonyms: 2'-fluoro-dAdo, NSC 274941
MF: $C_{10}H_{12}FN_5O_3$
FW: 269.2
Purity: $\geq 98\%$
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'-Fluoro-2'-deoxyadenosine is supplied as a solid. A stock solution may be made by dissolving the 2'-fluoro-2'-deoxyadenosine in the solvent of choice, which should be purged with an inert gas. 2'-Fluoro-2'-deoxyadenosine is soluble in DMSO. It is also soluble in water. We do not recommend storing the aqueous solution for more than one day.

Description

2'-Fluoro-2'-deoxyadenosine is a prodrug form of the cytotoxic purine nucleoside 2-fluoro-adenine.¹ *In vivo*, 2'-fluoro-2'-deoxyadenosine (100 mg/kg) reduces mean tumor weight in an *E. coli* purine nucleoside phosphorylase-expressing D54 glioblastoma mouse xenograft model.

Reference

1. Parker, W.B., Allan, P.W., Hassan, A.E., *et al.* Antitumor activity of 2-fluoro-2'-deoxyadenosine against tumors that express *Escherichia coli* purine nucleoside phosphorylase. *Cancer Gene Ther.* **10(1)**, 23-29 (2003).

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, 02/15/2024

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