

# PRODUCT INFORMATION

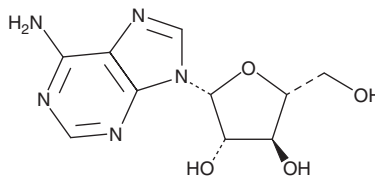


## Vidarabine

Item No. 18149

**CAS Registry No.:** 5536-17-4  
**Formal Name:** 9-β-D-arabinofuranosyl-9H-purin-6-amine  
**Synonyms:** Adenine Arabinoside, Ara-A, 9-β-D-Arabinofuranosyl adenine, Arabinosyladenine, NSC 247519, NSC 404241, Vira-A

**MF:** C<sub>10</sub>H<sub>13</sub>N<sub>5</sub>O<sub>4</sub>  
**FW:** 267.2  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 260 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

Vidarabine is supplied as a crystalline solid. A stock solution may be made by dissolving the vidarabine in the solvent of choice. Vidarabine is soluble in organic solvents such as DMSO and dimethyl formamide, which should be purged with an inert gas. The solubility of vidarabine in these solvents is approximately 11 and 3 mg/ml, respectively.

Vidarabine is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, vidarabine should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Vidarabine has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

Vidarabine is an analog of the nucleoside adenosine that has antiviral properties. It acts as a prodrug that, once phosphorylated by cellular enzymes, acts as both substrate and inhibitor of DNA polymerase.<sup>1</sup> Vidarabine is particularly effective against *H. simplex* and *V. zoster* viruses.<sup>1</sup>

### Reference

1. Bean, B. Antiviral therapy: Current concepts and practices. *Clin. Microbiol. Rev.* **5**(2), 146-182 (1992).

#### 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, 12/02/2022

#### 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](http://WWW.CAYMANCHEM.COM)