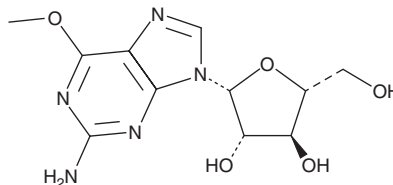


PRODUCT INFORMATION

Nelarabine

Item No. 20248

CAS Registry No.: 121032-29-9
Formal Name: 9-β-D-arabinofuranosyl-6-methoxy-9H-purin-2-amine
Synonyms: GW 506U78, Nelzarabine
MF: C₁₁H₁₅N₅O₅
FW: 297.3
Purity: ≥98%
UV/Vis.: λ_{max}: 211, 247, 281 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

Nelarabine is supplied as a crystalline solid. A stock solution may be made by dissolving the nelarabine in the solvent of choice. Nelarabine is soluble in organic solvents such as DMSO and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of nelarabine in these solvents is approximately 10 and 3 mg/ml, respectively.

Nelarabine is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, nelarabine should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Nelarabine 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

Nelarabine is a purine nucleoside analog that is used as a cytotoxic agent in hematological malignancies.¹ Nelarabine is converted into arabinosyl-guanine (ara-G) in the blood, and ara-G is selectively accumulated in T cells and malignant T-lymphoid cells, where it is phosphorylated to produce ara-GTP and incorporated into DNA.^{2,3} This leads to DNA breaks, increased expression of Fas ligand, and apoptosis.^{1,3} Ara-G is accumulated most effectively by acute lymphoid leukemia and chronic myeloid leukemia blasts.¹

References

1. Robak, T., Korycka, A., Lech-Maranda, E., *et al.* Current status of older and new purine nucleoside analogues in the treatment of lymphoproliferative diseases. *Molecules* **14**(3), 1183-1226 (2009).
2. Rodriguez, C.O., Jr., Mitchell, B.S., Ayres, M., *et al.* Arabinosylguanine is phosphorylated by both cytoplasmic deoxycytidine kinase and mitochondrial deoxyguanosine kinase. *Cancer Res.* **62**(11), 3100-3105 (2002).
3. Rodriguez, C.O., Jr., Stellrecht, C.M., and Gandhi, V. Mechanisms for T-cell selective cytotoxicity of arabinosylguanine. *Blood* **102**(5), 1842-1848 (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, 12/19/2022

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