

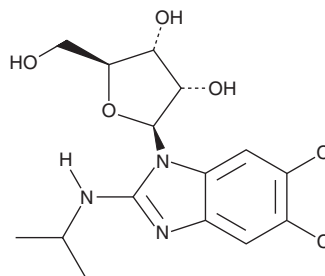
PRODUCT INFORMATION



Maribavir

Item No. 39650

CAS Registry No.: 176161-24-3
Formal Name: 5,6-dichloro-N-(1-methylethyl)-1-β-L-ribofuranosyl-1H-benzimidazol-2-amine
Synonyms: BW 1263W94, GW 257406X, MBV
MF: C₁₅H₁₉Cl₂N₃O₄
FW: 376.2
Purity: ≥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

Maribavir is supplied as a solid. A stock solution may be made by dissolving the maribavir in the solvent of choice, which should be purged with an inert gas. Maribavir is soluble in acetonitrile and DMSO.

Description

Maribavir is an antiviral agent.¹ It inhibits human cytomegalovirus (CMV) pUL97 protein kinase (IC₅₀ = 3 nM). Maribavir inhibits the replication of clinical isolates of CMV, including those resistant to ganciclovir (Item No. 37487), in infected MRC-5 human lung fibroblasts (IC_{50s} = 0.03-0.13 μM). Oral administration of maribavir (25 or 75 mg/kg) decreases CMV titers in a mouse model of ocular CMV infection following implantation of human fetal retinal tissue.² Formulations containing maribavir have been used in the treatment of post-transplant CMV infections.

References

1. Biron, K.K., Harvey, R.J., Chamberlain, S.C., *et al.* Potent and selective inhibition of human cytomegalovirus replication by 1263W94, a benzimidazole L-riboside with a unique mode of action. *Antimicrob. Agents Chemother.* **46(8)**, 2365-2372 (2002).
2. Kern, E.R., Hartline, C.B., Rybak, R.J., *et al.* Activities of benzimidazole D- and L-ribonucleosides in animal models of cytomegalovirus infections. *Antimicrob. Agents Chemother.* **48(5)**, 1749-1755 (2004).

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, 01/30/2024

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