

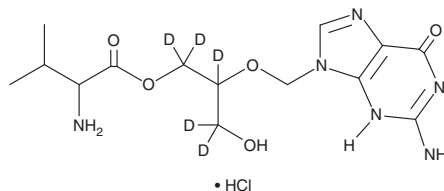
PRODUCT INFORMATION



Valganciclovir-d₅ (hydrochloride)

Item No. 30746

Formal Name: DL-valine, 2-[(2-amino-1,6-dihydro-6-oxo-9H-purin-9-yl)methoxy]-3-hydroxypropyl-1,1,2,3,3-d₅ ester, monohydrochloride
MF: C₁₄H₁₇D₅N₆O₅ • HCl
FW: 395.9
Chemical Purity: ≥95% (Valganciclovir)
Deuterium Incorporation: ≥99% deuterated forms (d₁-d₅); ≤1% d₀
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

Valganciclovir-d₅ (hydrochloride) is intended for use as an internal standard for the quantification of valganciclovir (Item No. 13875) by GC- or LC-MS. The accuracy of the sample weight in this vial is between 5% over and 2% under the amount shown on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard by constructing a standard curve of peak intensity ratios (deuterated versus unlabeled).

Valganciclovir-d₅ (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the valganciclovir-d₅ (hydrochloride) in the solvent of choice, which should be purged with an inert gas. Valganciclovir-d₅ (hydrochloride) is soluble in organic solvents such as methanol and DMSO.

Description

Valganciclovir is a valyl ester prodrug form of the antiviral nucleoside analog ganciclovir (Item No. 13853). *In vivo*, valganciclovir (200 mg/kg) increases survival and reduces weight loss in an immunosuppressed hamster model of disseminated adenovirus type 5 (Ad5) infection.¹ Formulations containing valganciclovir have been used in the prevention of post-transplant cytomegalovirus (CMV) infections.

Reference

1. Toth, K., Ying, B., Tollefson, A.E., *et al.* Valganciclovir inhibits human adenovirus replication and pathology in permissive immunosuppressed female and male Syrian hamsters. *Viruses* **7**(3), 1409-1428 (2015).

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, 12/09/2022

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