

# PRODUCT INFORMATION



## Valganciclovir (hydrochloride)

Item No. 13875

**CAS Registry No.:** 175865-59-5  
**Formal Name:** 2-[(2-amino-1,6-dihydro-6-oxo-9H-purin-9-yl)methoxy]-3-hydroxypropyl ester L-valine, monohydrochloride

**Synonym:** Ro 107-9070/194

**MF:** C<sub>14</sub>H<sub>22</sub>N<sub>6</sub>O<sub>5</sub> • HCl

**FW:** 390.8

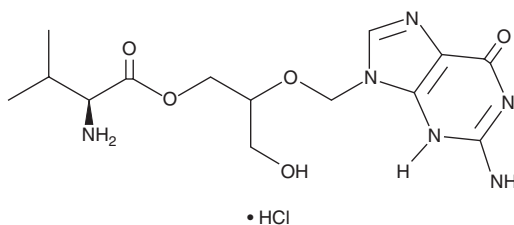
**Purity:** ≥95%

**UV/Vis.:** λ<sub>max</sub>: 255 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

Valganciclovir (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the valganciclovir (hydrochloride) in the solvent of choice, which should be purged with an inert gas. Valganciclovir (hydrochloride) is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of valganciclovir (hydrochloride) in these solvents is approximately 20 and 1 mg/ml, respectively.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of valganciclovir (hydrochloride) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of valganciclovir (hydrochloride) in PBS (pH 7.2) is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### 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.<sup>1</sup> 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, 11/30/2022

#### 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