

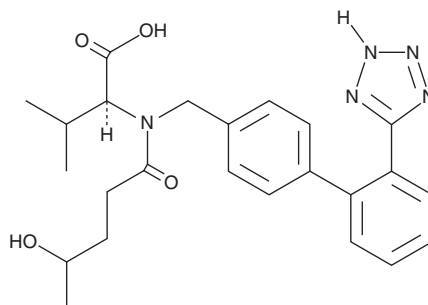
# PRODUCT INFORMATION



## 4-hydroxy Valsartan

Item No. 27849

**CAS Registry No.:** 188259-69-0  
**Formal Name:** N-(4-hydroxy-1-oxopentyl)-N-[[2'-(2H-tetrazol-5-yl)[1,1'-biphenyl]-4-yl]methyl]-L-valine  
**Synonyms:** CGP 71580, Valeryl 4-hydroxy Valsartan  
**MF:** C<sub>24</sub>H<sub>29</sub>N<sub>5</sub>O<sub>4</sub>  
**FW:** 451.5  
**Purity:** ≥98% (mixture of diastereomers)  
**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

4-hydroxy Valsartan is supplied as a solid. A stock solution may be made by dissolving the 4-hydroxy valsartan in the solvent of choice, which should be purged with an inert gas. 4-hydroxy Valsartan is slightly soluble in methanol.

### Description

4-hydroxy Valsartan is a major metabolite of the angiotensin II type 1 (AT<sub>1</sub>) receptor antagonist valsartan (Item No. 14178).<sup>1</sup> It reduces platelet aggregation induced by epinephrine (Item No. 21245) and collagen but not ADP in human whole blood.

### References

1. Serebruany, V.L., Malinin, A.I., Lowry, D.R., *et al.* Effects of valsartan and valeryl 4-hydroxy valsartan on human platelets: A possible additional mechanism for clinical benefits. *J. Cardiovasc. Pharmacol.* **43**(5), 677-684 (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, 10/14/2022

#### 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