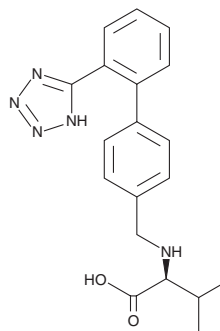


PRODUCT INFORMATION

Desvaleryl Valsartan

Item No. 35634

CAS Registry No.: 676129-92-3
Formal Name: N-[[2'-(2H-tetrazol-5-yl)[1,1'-biphenyl]-4-yl]methyl]-L-valine
Synonym: Valsartan Impurity B
MF: C₁₉H₂₁N₅O₂
FW: 351.4
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

Desvaleryl valsartan is supplied as a solid. A stock solution may be made by dissolving the desvaleryl valsartan in the solvent of choice, which should be purged with an inert gas. Desvaleryl valsartan is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of desvaleryl valsartan in these solvents is approximately 1, 5, and 3 mg/ml, respectively.

Description

Desvaleryl valsartan is a potential impurity found in commercial preparations of the angiotensin II type 1 (AT₁) receptor antagonist valsartan.¹ It is a degradation product formed under acidic conditions.²

References

1. Krishnaiah, C., Reddy, A.R., Kumar, R., *et al.* Stability-indicating UPLC method for determination of valsartan and their degradation products in active pharmaceutical ingredient and pharmaceutical dosage forms. *J. Pharm. Biomed. Anal.* **53**(3), 483-489 (2010).
2. Kumaraswamy, K., Gandham, H.B., Prasad, R.J.D., *et al.* Isolation and characterization of novel degradation products of valsartan by NMR and high resolution mass spectroscopy: Development and validation of valsartan by UPLC. *Asian J. Chem.* **32**(5), 1064-1068 (2020).

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, 11/14/2022

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