

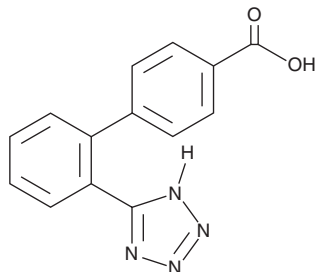
PRODUCT INFORMATION



Valsartan Acid

Item No. 35361

CAS Registry No.: 164265-78-5
Formal Name: 2'-(2H-tetrazol-5-yl)-[1,1'-biphenyl]-4-carboxylic acid
MF: $C_{14}H_{10}N_4O_2$
FW: 266.3
Purity: $\geq 95\%$
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

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

Description

Valsartan acid is a transformation product of the angiotensin II type 1 (AT_1) receptor antagonist valsartan (Item No. 14178).¹ It is formed during activated sludge treatment and has been found as a contaminant in drinking water.

Reference

1. Nödler, K., Hillebrand, O., Idzik, K., *et al.* Occurrence and fate of the angiotensin II receptor antagonist transformation product valsartan acid in the water cycle--a comparative study with selected β -blockers and the persistent anthropogenic wastewater indicators carbamazepine and acesulfame. *Water Res.* **47**(17), 6650-6659 (2013).

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, 09/30/2022

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