

PRODUCT INFORMATION



5-(4'-(Azidomethyl)(1,1'-biphenyl)-2-yl)-2H-tetrazole

Item No. 38780

CAS Registry No.: 152708-24-2

Synonym: AZTT

MF: $C_{14}H_{11}N_7$

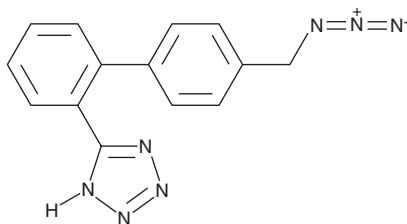
FW: 277.3

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

5-(4'-(Azidomethyl)(1,1'-biphenyl)-2-yl)-2H-tetrazole (AZTT) is supplied as a solid. A stock solution may be made by dissolving the AZTT in the solvent of choice, which should be purged with an inert gas. AZTT is soluble in methanol and DMSO.

Description

AZTT is a potential impurity found in commercial preparations of tetrazole-containing sartans, including losartan, irbesartan, and valsartan.¹

Reference

1. Wichitnithad, W., Nantaphol, S., Thitikornpong, W., *et al.* Development and validation of an LC-MS/MS method for simultaneous determination of three organic azido impurities in tetrazole-containing sartans. *Arab. J. Chem.* **16**(8), 104951 (2023).

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, 12/08/2023

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