

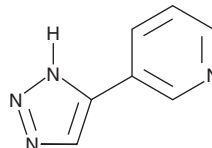
PRODUCT INFORMATION



3-TYP

Item No. 29660

CAS Registry No.: 120241-79-4
Formal Name: 3-(1H-1,2,3-triazol-5-yl)-pyridine
MF: C₇H₆N₄
FW: 146.2
Purity: ≥98%
UV/Vis.: λ_{max}: 245 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

3-TYP is supplied as a crystalline solid. A stock solution may be made by dissolving the 3-TYP in the solvent of choice, which should be purged with an inert gas. 3-TYP is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of 3-TYP in ethanol is approximately 5 mg/ml and approximately 10 mg/ml in DMSO and DMF.

3-TYP is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, 3-TYP should first be dissolved in DMF and then diluted with the aqueous buffer of choice. 3-TYP has a solubility of approximately 0.16 mg/ml in a 1:5 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

3-TYP is a sirtuin 3 (SIRT3) inhibitor (IC₅₀ = 38 μM).¹ It is selective for SIRT3 over SIRT1 and SIRT2 at 1 mM. 3-TYP (1 mM) increases acetylation of lysine residues in mitochondrial proteins in HeLa cells. It inhibits the growth of CCRF-CEM, COLO 205, OVCAR-4, OVCAR-8, and SKOV3 cells by greater than 50% when used at a concentration of 10 μM.

Reference

1. Galli, U., Mesenzani, O., Coppo, C., *et al.* Identification of a sirtuin 3 inhibitor that displays selectivity over sirtuin 1 and 2. *Eur. J. Med. Chem.* **55**, 58-66 (2012).

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, 12/09/2022

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