

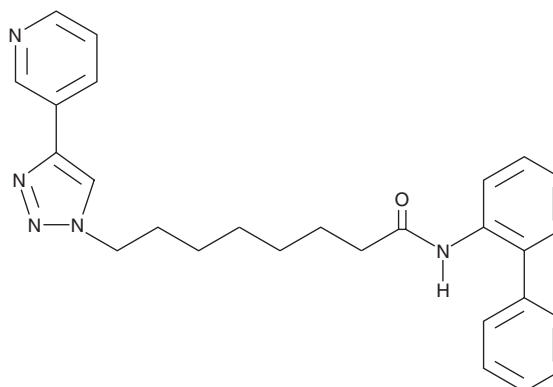
PRODUCT INFORMATION



GPP 78

Item No. 13670

CAS Registry No.: 1202580-59-3
Formal Name: N-[1,1'-biphenyl]-2-yl-4-(3-pyridinyl)-1H-1,2,3-triazole-1-octanamide
Synonyms: CAY10618, NMPRTase Inhibitor
MF: C₂₇H₂₉N₅O
FW: 439.6
Purity: ≥95%
UV/Vis.: λ_{max}: 234 nm
Supplied as: A solution in methanol
Storage: -20°C
Stability: ≥2 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

GPP 78 is supplied as a solution in methanol. To change the solvent, simply evaporate the GPP 78 under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as ethanol, DMSO, and dimethyl formamide purged with an inert gas can be used. The solubility of GPP 78 in these solvents is approximately 10 mg/ml.

GPP 78 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the ethanolic solution of GPP 78 should be diluted with the aqueous buffer of choice. GPP 78 has a solubility of approximately 0.25 mg/ml in a 1:3 solution of DMSO:PBS (pH 7.2) using this method.

Description

GPP 78 is a potent inhibitor of Nampt (IC₅₀ = 3 nM).¹ It induces cell death in the neuroblastoma cell line SH-SY5Y with an IC₅₀ value of 3.8 nM through a process that appears to involve autophagy.²

References

- Colombano, G., Travelli, C., Galli, U., *et al.* A novel potent nicotinamide phosphoribosyltransferase inhibitor synthesized via click chemistry. *J. Med. Chem.* **53**(2), 616-623 (2010).
- Stein, R.L., Melandri, F., and Dick, L. Kinetic characterization of the chymotryptic activity of the 20S proteasome. *Biochemistry* **35**(13), 3899-3908 (1996).

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, 07/24/2018

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