

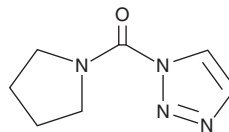
PRODUCT INFORMATION



AA26-8

Item No. 36937

CAS Registry No.: 1312782-34-5
Formal Name: 1-pyrrolidinyl-1H-1,2,3-triazol-1-yl-methanone
Synonyms: AA26-9, KK007-N1
MF: C₇H₁₀N₄O
FW: 166.2
Purity: ≥98%
UV/Vis.: λ_{max}: 218 nm
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

AA26-8 is supplied as a solid. A stock solution may be made by dissolving the AA26-8 in the solvent of choice, which should be purged with an inert gas. AA26-8 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of AA26-8 in these solvents is approximately 3, 5, and 10 mg/ml, respectively.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of AA26-8 can be prepared by directly dissolving the solid in aqueous buffers. The solubility of AA26-8 in PBS (pH 7.2) is approximately 2 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

AA26-8 is a pan-serine hydrolase inhibitor.¹ It inhibits acylaminoacyl-peptide hydrolase (APEH), platelet-activating factor acetylhydrolase 2 (PAF-AH2), α/β hydrolase domain-containing protein 11 (ABHD11), lysophospholipase 1 (LYPLA1), fatty acid amide hydrolase (FAAH), and neutral cholesterol ester hydrolase 1 (NCEH1) when used at a concentration of 1 μM. AA26-8 (10 μM) also inhibits the interaction of the α/β hydrolase Dwarf14 (D14) and a negative regulator of its signaling, D53, induced by the plant growth regulator GR24 (Item No. 13210) in a yeast two-hybrid assay.²

References

1. Adibekian, A., Martin, B.R., Wang, C., *et al.* Click-generated triazole ureas as ultrapotent *in vivo*-active serine hydrolase inhibitors. *Nat. Chem. Biol.* **7**(7), 469-478 (2011).
2. Nakamura, H., Hirabayashi, K., Miyakawa, T., *et al.* Triazole ureas covalently bind to strigolactone receptor and antagonize strigolactone responses. *Mol. Plant* **12**(1), 44-58 (2019).

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, 12/21/2022

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