

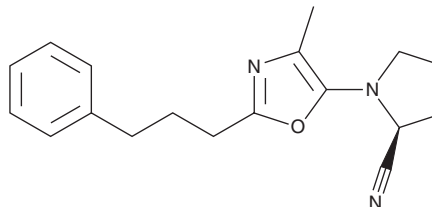
PRODUCT INFORMATION



HUP-55

Item No. 39117

CAS Registry No.: 3006795-68-9
Formal Name: (2S)-1-[4-methyl-2-(3-phenylpropyl)-5-oxazolyl]-2-pyrrolidinecarbonitrile
MF: C₁₈H₂₁N₃O
FW: 295.4
Purity: ≥98%
Supplied as: A 10 mg/ml solution in acetonitrile
Storage: -20°C
Stability: ≥2 years
Special Conditions: Oxygen and light-sensitive



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

Description

HUP-55 is an inhibitor of prolyl endopeptidase (IC₅₀ = 5 nM).¹ It reduces α-synuclein dimerization in Neuro2a cells, induces autophagy in HEK293 cells, and decreases hydrogen peroxide-induced increases in reactive oxygen species (ROS) production in SH-SY5Y cells when used at a concentration of 10 μM. HUP-55 (10 mg/kg) decreases ipsilateral paw use and striatal oligomeric α-synuclein levels in a mouse model of Parkinson's disease induced by adenoviral delivery of an α-synuclein-encoding vector (AAV-αSyn).

Reference

1. Kilpeläinen, T.P., Päätsi, H.T., Svarcbahts, R., *et al.* Nonpeptidic oxazole-based prolyl oligopeptidase ligands with disease-modifying effects on α-synuclein mouse models of Parkinson's disease. *J. Med. Chem.* **66**(11), 7475-7496 (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, 01/26/2026

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