

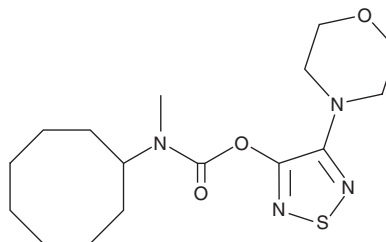
PRODUCT INFORMATION



JZP 430

Item No. 35383

CAS Registry No.: 1672691-74-5
Formal Name: N-cyclooctyl-N-methyl-carbamic acid, 4-(4-morpholinyl)-1,2,5-thiadiazol-3-yl ester
MF: C₁₆H₂₆N₄O₃S
FW: 354.5
Purity: ≥98%
UV/Vis.: λ_{max}: 206, 210, 302 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

JZP 430 is supplied as a solid. A stock solution may be made by dissolving the JZP 430 in the solvent of choice, which should be purged with an inert gas. JZP 430 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of JZP 430 in DMF is approximately 30 mg/ml and approximately 10 mg/ml in ethanol and DMSO.

Description

JZP 430 is an inhibitor of α/β-hydrolase domain-containing protein 6 (ABHD6; IC₅₀ = 0.044 for the human enzyme).¹ It is selective for ABHD6 over fatty acid amid hydrolase (FAAH) and monoacylglycerol lipase (MAGL) when used at a concentration of 2.5 μM, ABHD12 at 1 μM, as well as lysosomal acid lipase (LAL) and cannabinoid 1 (CB₁) and CB₂ receptors at 10 μM.

Reference

1. Patel, J.Z., Nevalainen, T.J., Savinainen, J.R., *et al.* Optimization of 1,2,5-thiadiazole carbamates as potent and selective ABHD6 inhibitors. *ChemMedChem* **10**(2), 253-265 (2015).

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, 11/03/2022

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