

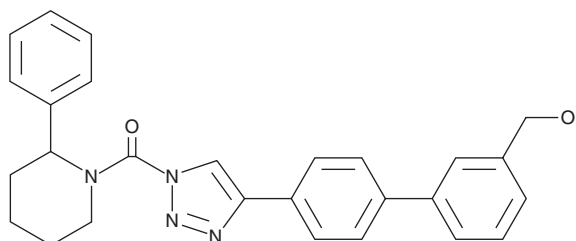
PRODUCT INFORMATION



KT182

Item No. 15404

CAS Registry No.: 1402612-62-7
Formal Name: [4-[3'-(hydroxymethyl)[1,1'-biphenyl]-4-yl]-1H-1,2,3-triazol-1-yl](2-phenyl-1-piperidiny)-methanone
MF: C₂₇H₂₆N₄O₂
FW: 438.5
Purity: ≥98%
UV/Vis.: λ_{max}: 280 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

KT182 is supplied as a crystalline solid. A stock solution may be made by dissolving the KT182 in the solvent of choice. KT182 is soluble in organic solvents such as DMSO and dimethyl formamide, which should be purged with an inert gas. The solubility of KT182 in these solvents is approximately 10 mg/ml.

Description

KT182 is a potent inhibitor of α/β -hydrolase domain-containing protein 6 (ABHD6) with IC₅₀ values of 1.7, 15.1, and 0.24 nM using Neuro2a membrane proteomes, recombinant ABHD6 in HEK293T cells, and Neuro2a cells *in situ*, respectively.¹ Following administration of KT182 (1 mg/kg, i.p.) in mice, ABHD6 is inactivated in liver and brain extracts, suggesting that it is brain-penetrant, unlike the closely related compound KT203 (Item No. 14819).

Reference

1. Hsu, K.L., Tsuboi, K., Chang, J.W., *et al.* Discovery and optimization of piperidyl-1,2,3-triazole ureas as potent, selective, and *in vivo*-active inhibitors of α/β -hydrolase domain containing 6 (ABHD6). *J. Med. Chem.* **56**(21), 8270-8279 (2013).

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, 09/11/2017

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