

# PRODUCT INFORMATION



## KT203

Item No. 14819

CAS Registry No.: 1402612-64-9

Formal Name: 4'-[1-[[2-(phenylmethyl)-1-piperidinyl]carbonyl]-1H-1,2,3-triazol-4-yl]-[1,1'-biphenyl]-3-carboxylic acid

Synonym: ML-296

MF: C<sub>28</sub>H<sub>26</sub>N<sub>4</sub>O<sub>3</sub>

FW: 466.5

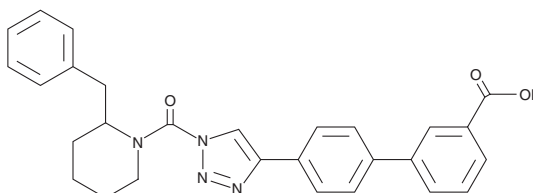
Purity: ≥95%

UV/Vis.: λ<sub>max</sub>: 277 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

KT203 is supplied as a crystalline solid. A stock solution may be made by dissolving the KT203 in the solvent of choice. KT203 is soluble in organic solvents such as DMSO and dimethyl formamide, which should be purged with an inert gas. The solubility of KT203 in these solvents is approximately 10 mg/ml.

### Description

KT203 is a potent inhibitor of  $\alpha/\beta$ -hydrolase domain-containing protein 6 (ABHD6) with IC<sub>50</sub> values of 0.82, 3.9, and 0.31 nM using Neuro2A membrane proteomes, recombinant ABHD6 in HEK293T cells, and Neuro2A cells *in situ*, respectively.<sup>1</sup> Following administration of KT203 (1 mg/kg, i.p.) in mice, ABHD6 is inactivated in liver but not brain extracts, suggesting that KT203 is not brain-penetrant.

### 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  $\alpha/\beta$ -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, 08/15/2017

#### 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