

PRODUCT INFORMATION

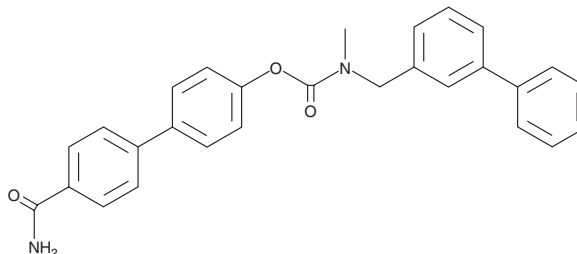


WWL123

Item No. 16849

CAS Registry No.: 1338574-83-6
Formal Name: N-([1,1'-biphenyl]-3-ylmethyl)-N-methyl-carbamic acid, 4'-(aminocarbonyl) [1,1'-biphenyl]-4-yl ester

MF: C₂₈H₂₄N₂O₃
FW: 436.5
Purity: ≥95%
UV/Vis.: λ_{max}: 265 nm
Supplied as: A crystalline 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

WWL123 is supplied as a crystalline solid. A stock solution may be made by dissolving the WWL123 in the solvent of choice. WWL123 is soluble in the organic solvent DMSO, which should be purged with an inert gas, at a concentration of approximately 1 mg/ml.

WWL123 is sparingly soluble in aqueous solutions. To enhance aqueous solubility, dilute the organic solvent solution into aqueous buffers or isotonic saline. If performing biological experiments, ensure the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

Description

The serine hydrolase known as α/β-hydrolase domain-containing protein 6 (ABHD6) hydrolyzes 2-arachidonoyl glycerol (Item No. 62160) to regulate its availability at cannabinoid receptors.¹ WWL123 is a brain-penetrant inhibitor of ABHD6 (IC₅₀ = 0.43 μM) that demonstrates >10-fold selectivity for ABHD6 compared to a panel of ~35 other serine hydrolases.² Inhibition of ABHD6 by WWL123 has been used to decrease seizure incidence in a genetic mouse model of juvenile Huntington's disease as well as in chemically-induced epilepsy models.³

References

- Blankman, J.L., Simon, G.M., and Cravatt, B.F. A comprehensive profile of brain enzymes that hydrolyze the endocannabinoid 2-arachidonoylglycerol. *Chem. Biol.* **14**, 1347-1356 (2007).
- Bachovchin, D.A., Ji, T., Simon, G.M., et al. Superfamily-wide portrait of serine hydrolase inhibition achieved by library-versus-library screening. *Proc. Natl. Acad. Sci. USA* **107**(49), 20941-20946 (2010).
- Naydenov, A.V., Horne, E.A., Cheah, C.S., et al. ABHD6 blockade exerts antiepileptic activity in PTZ-induced seizures and in spontaneous seizures in R6/2 mice. *Neuron* **83**(2), 361-371 (2014).

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, 02/09/2024

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