

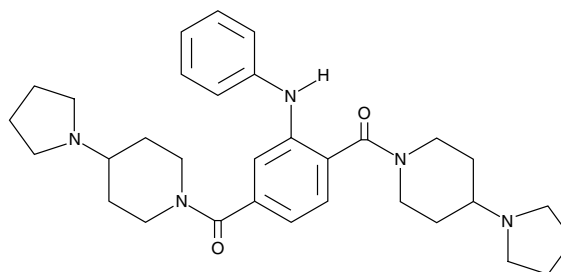
Product Information



UNC1215

Item No. 13968

CAS Registry No.: 1415800-43-9
Formal Name: (2-(phenylamino)-1,4-phenylene)bis((4-(pyrrolidin-1-yl)piperidin-1-yl)methanone)
MF: C₃₂H₄₃N₅O₂
FW: 529.7
Purity: ≥98%
Stability: ≥2 years at -20°C
Supplied as: A crystalline solid
UV/Vis.: λ_{max}: 289 nm



Laboratory Procedures

For long term storage, we suggest that UNC1215 be stored as supplied at -20°C. It should be stable for at least two years.

UNC1215 is supplied as a crystalline solid. A stock solution may be made by dissolving the UNC1215 in the solvent of choice. UNC1215 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of UNC1215 in these solvents is approximately 14, 5, and 16 mg/ml, respectively.

UNC1215 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, UNC1215 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. UNC1215 has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Methyllysine (Kme) recognition “reader” domains play a central role in epigenetic regulation during cellular differentiation, development, and gene transcription. UNC1215 is a potent and selective chemical probe for the Kme reading function of L3MBTL3, a member of the malignant brain tumor (MBT) family of chromatin interacting transcriptional repressors.¹ UNC1215 binds L3MBTL3 with a K_d value of 120 nM (IC₅₀ = 40 nM), competitively displacing mono- or dimethyl-lysine containing peptides.¹ This probe is greater than 50-fold selective toward L3MBTL3 than other members of the human MBT family and demonstrates selectivity against more than 200 other Kme reader domains examined.¹

Reference

1. James, L.I., Barsyte-Lovejoy, D., Zhong, N., *et al.* Discovery of a chemical probe for the L3MBTL3 methyllysine reader domain. *Nat. Chem. Biol.* [In press] (2013).

Related Products

For a list of related products please visit: www.caymanchem.com/catalog/13968

WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY: NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

MATERIAL SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. Before use, the user must review the complete Material Safety Data Sheet, which has been sent *via* email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman **warrants only** to the original customer that the material will **meet our specifications at the time of delivery.**

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have **any obligation or liability**, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees.

Buyer's **exclusive remedy** and Cayman's sole liability hereunder shall be limited to a **refund** of the purchase price, or at Cayman's option, the **replacement**, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our **Warranty and Limitation of Remedy located on our website and in our catalog.**

Copyright Cayman Chemical Company, 01/30/2013

Cayman Chemical

Mailing address

1180 E. Ellsworth Road
Ann Arbor, MI
48108 USA

Phone

(800) 364-9897
(734) 971-3335

Fax

(734) 971-3640

E-Mail

custserv@caymanchem.com

Web

www.caymanchem.com