

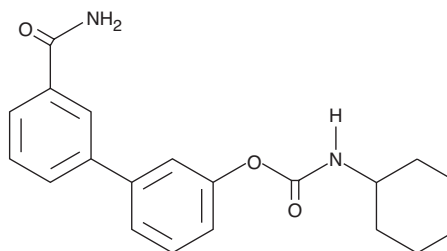
Product Information



URB597

Catalog No. 10046

CAS Registry No.: 546141-08-6
Formal Name: (3'-(aminocarbonyl)[1,1'-biphenyl]-3-yl)-cyclohexylcarbamate
MF: C₂₀H₂₂N₂O₃
FW: 338.4
Purity: ≥98%
Stability: ≥2 years at -20°C
Supplied as: A crystalline solid



Laboratory Procedures

For long term storage, we suggest that URB597 be stored as supplied at -20°C. It will be stable for at least two years.

URB597 is supplied as a crystalline solid. A stock solution may be made by dissolving the URB597 in an organic solvent purged with an inert gas. URB597 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of URB597 in ethanol is approximately 1 mg/ml, and 10 mg/ml in DMSO and DMF.

URB597 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, URB597 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. URB597 has a solubility of 500 µg/ml in a 1:1 solution of DMF:PBS (pH 7.2) or in a 1:2 solution of DMSO:PBS (pH 7.2) using this method. Therefore, further dilutions of the organic solvent solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that 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.

The enzyme fatty acid amide hydrolase (FAAH) is widely expressed in brain and other tissues, and is capable of hydrolyzing anandamide (AEA) and other simple esters and amides with long unsaturated acyl chains.¹ URB597 was recently synthesized and reported to be a potent and selective inhibitor of FAAH.² URB597 inhibits FAAH with an IC₅₀ of 4.6 nM in brain membranes and 0.5 nM in intact neurons. It has been demonstrated in FAAH^(-/-) mice that marked depression of FAAH activity results in reduced sensation of pain and enhanced endocannabinoid signalling.³ URB597 exhibits both anti-nociceptive and anxiolytic effects *in vivo* without evoking other symptoms associated with cannabinoid-like compounds. Thus, URB597 may serve as a lead compound for the development of new analgesic and anxiolytic drugs.

References

1. Cravatt, B.F., Giang, D.K., Mayfield, S.P., *et al.* Molecular characterization of an enzyme that degrades neuromodulatory fatty-acid amides. *Nature* **384**, 83-87 (1996).
2. Kathuria, S., Gaetani, S., Fegley, D., *et al.* Modulation of anxiety through blockade of anandamide hydrolysis. *Nature Med.* **1(9)**, 76-81 (2003).
3. Cravatt, B.F., Demarest, K., Patricelli, M.P., *et al.* Supersensitivity to anandamide and enhanced endogenous cannabinoid signaling in mice lacking fatty acid amide hydrolase. *Proc. Natl. Acad. Sci. USA* **98(16)**, 9371-9376 (2001).

Related Products

Methyl α-Linolenyl Fluorophosphonate - Cat. No. 70662 • Methyl γ-Linolenyl Fluorophosphonate - Cat. No. 70664 • Methyl Nonadecadienyl Fluorophosphonate - Cat. No. 70665 • CAY10400 - Cat. No. 71650

WARNING: THIS PRODUCT IS NOT INTENDED OR APPROVED FOR HUMAN OR VETERINARY USE. USE OF THIS PRODUCT FOR HUMAN OR ANIMAL TESTING IS EXTREMELY HAZARDOUS AND MAY RESULT IN DISEASE, SEVERE INJURY, OR DEATH.

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 under separate cover to the MSDS supervisor at 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, 09/08/2006

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