

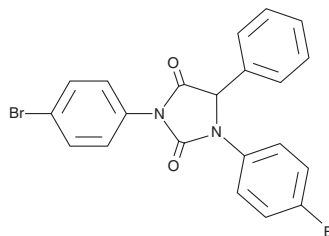
Product Information



CAY10508

Item No. 10008669

CAS Registry No.: 878533-35-8
Formal Name: 1,3-bis(4-bromophenyl)-5-phenyl-2,4-imidazolidinedione
MF: C₂₁H₁₄Br₂N₂O₂
FW: 486.2
Purity: ≥98%
Stability: ≥2 years at -20°C
Supplied as: A crystalline solid



Laboratory Procedures

For long term storage, we suggest that CAY10508 be stored as supplied at -20°C. It should be stable for at least two years.

CAY10508 is supplied as a crystalline solid. A stock solution may be made by dissolving the CAY10508 in an organic solvent purged with an inert gas. CAY10508 is soluble in organic solvents such as DMSO and dimethyl formamide (DMF). The solubility of CAY10508 in DMSO is approximately 10 mg/ml and approximately 30 mg/ml in DMF.

CAY10508 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, CAY10508 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. CAY10508 has a solubility of approximately 0.25 mg/ml in a 1:2 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Central cannabinoid (CB₁) receptor antagonists and inverse agonists exhibit therapeutic potential for obesity and drug dependence, while being devoid of psychotropic effects. CAY10508 is a potent and selective CB₁ receptor inverse agonist with a K_i value of 243 nM and an EC₅₀ of 195 nM.¹ A 10 μM concentration of CAY10508 resulted in 100% and 35% displacement of [³H]-CP-55,940 at the CB₁ and peripheral cannabinoid (CB₂) receptors, respectively. A [35S]-GTP_γS binding assay revealed the inverse agonist properties of CAY10508 at the CB₁ receptor.

Reference

1. Muccioli, G.G., Wouters, J., Charlier, C., *et al.* Synthesis and activity of 1,3,5-triphenylimidazolidine-2,4-diones and 1,3,5-triphenyl-2-thioxoimidazolidin-4-ones: Characterization of new CB₁ cannabinoid receptor inverse agonists/antagonists. *J. Med. Chem.* **49**, 872-882 (2006).

Related Products

LH 21 - Item No. 13453 • 2-Arachidonoyl Glycerol - Item No. 62160 • AM251 - Item No. 71670 • Arachidonoyl Ethanolamide - Item No. 90050 • O-Arachidonoyl Ethanolamine (hydrochloride) - Item No. 91050 • CB₁ Receptor (C-Term) Polyclonal Antibody - Item No. 10006590

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, 03/31/2011

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