

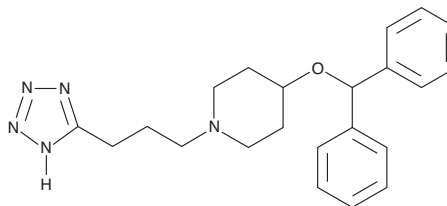
Product Information



HQL-79

Catalog No. 10134

CAS Registry No.: 162641-16-9
Formal Name: 4-(diphenylmethoxy)-1-[3-(1H-tetrazol-5-yl)propyl]-piperidine
MF: C₂₂H₂₇N₅O
FW: 377.5
Purity: ≥98%
Stability: ≥1 year at -20°C
Supplied as: A crystalline solid



Laboratory Procedures

For long term storage, we suggest that HQL-79 be stored as supplied at -20°C. It will be stable for at least one year.

HQL-79 is supplied as a crystalline solid. A stock solution may be made by dissolving the HQL-79 in an organic solvent purged with an inert gas. HQL-79 is soluble in organic solvents such as methanol. The solubility of HQL-79 in methanol is approximately 2.5 mg/ml.

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.

HQL-79 is a selective inhibitor of hematopoietic prostaglandin D (PGD) synthase. Structurally, it is a synthetic tetrazole compound originally prepared as a possible antihistamine.¹ Subsequent evaluation in models of allergy and asthma demonstrates the HQL-79 is orally available, and inhibits the synthesis of PGD₂.² HQL-79 is thus a likely lead compound for the preparation of potent, selective hematopoietic PGD synthase inhibitors.

References

1. Matsushita, K., Nobutoshi, A., and Aritake, K. Pharmacological studies on the novel antiallergic drug HQL-79: II. Elucidation of mechanisms for antiallergic and antiasthmatic effects. *Jpn. J. Pharmacol.* **78(1)**, 11-22 (1998).
2. Matsushita, K., Nobutoshi, A., and Hizue, A. Pharmacological studies on the novel antiallergic drug HQL-79: I. Antiallergic and antiasthmatic effects in various experimental models. *Jpn. J. Pharmacol.* **78(1)**, 1-10 (1998).

Related Products

Prostaglandin D₂ - Cat. No. 12010 • Prostaglandin D Synthase (Lipocalin-type) Polyclonal Antibody - Cat. No. 160003 • Prostaglandin D Synthase (hematopoietic) Polyclonal Antibody - Cat. No. 160013 • Prostaglandin E₂ Express EIA Kit - Cat. No. 500141 • Prostaglandin D₂-MOX Express EIA Kit - Cat. No. 500151 • Prostaglandin D₂-MOX EIA Kit - Cat. No. 512011

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 MSD supervisor at your institution.

WARRANTY AND LIMITATION OF REMEDY

Cayman Chemical Company makes no warranty of any kind, expressed or implied, including, but not limited to, the warranties of fitness for a particular purpose and merchantability, which extends beyond the description of the chemicals on the face hereof, except that the material will meet our specifications at the time of delivery. Buyer's exclusive remedy and Cayman Chemical Company's sole liability hereunder shall be limited to refund of the purchase price of, or at Cayman Chemical Company's option the replacement of, all material that does not meet our specifications. Cayman Chemical Company shall not be liable otherwise or for incidental or consequential damages, including, but not limited to, the costs of handling. Said refund or replacement is conditioned on Buyer giving written notice to Cayman Chemical Company within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within said thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

Copyright Cayman Chemical Company, 03/20/2006

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