

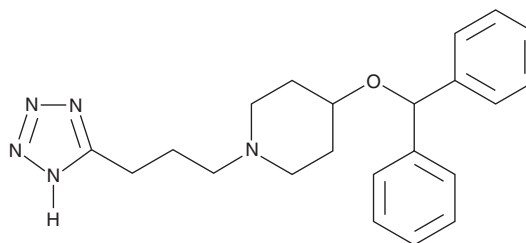
PRODUCT INFORMATION



HQL-79

Item 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%
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

HQL-79 is supplied as a crystalline solid. A stock solution may be made by dissolving the HQL-79 in the solvent of choice, which should be purged with an inert gas. HQL-79 is soluble in the organic solvent methanol at a concentration of approximately 2.5 mg/ml.

Description

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 (Item No. 12010).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).

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, 04/15/2024

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