

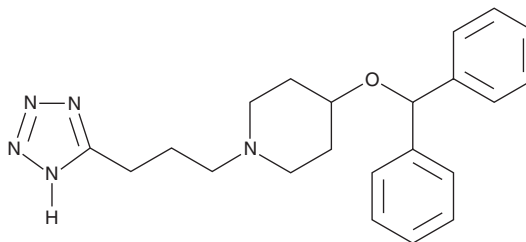
# PRODUCT INFORMATION



## HQL-79

Item No. 10134

CAS Registry No.: 162641-16-9  
Formal Name: 4-(diphenylmethoxy)-1-[3-(1H-tetrazol-5-yl)propyl]-piperidine  
MF: C<sub>22</sub>H<sub>27</sub>N<sub>5</sub>O  
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 should 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.

### 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.<sup>1</sup> Subsequent evaluation in models of allergy and asthma demonstrates the HQL-79 is orally available, and inhibits the synthesis of PGD<sub>2</sub> (Item No. 12010).<sup>2</sup> 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, 08/13/2015

#### 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