

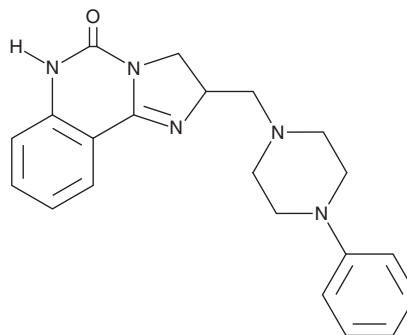
PRODUCT INFORMATION



2-PMDQ

Item No. 43206

CAS Registry No.: 139047-55-5
Formal Name: 2,6-dihydro-2-[(4-phenyl-1-piperazinyl)methyl]-imidazo[1,2-c]quinazolin-5(3H)-one
Synonym: AT-112
MF: C₂₁H₂₃N₅O
FW: 361.4
Purity: ≥98%
Supplied as: A 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

2-PMDQ is supplied as a solid. A stock solution may be made by dissolving the 2-PMDQ in the solvent of choice, which should be purged with an inert gas. 2-PMDQ is slightly soluble (0.1-1 mg/ml) in DMSO.

Description

2-PMDQ is an antagonist of α_1 -adrenergic receptors (α_1 -ARs; $K_i = 1.7$ nM).¹ It is selective for α_1 -ARs over α_2 -ARs ($K_i = 962$ nM). 2-PMDQ has antihypertensive activity in anesthetized rats ($ED_{50} = 0.25$ mg/kg).²

References

1. Chern, J.W., Tao, P.L., Yen, M.H., *et al.* Studies on quinazolines. 5. 2,3-dihydroimidazo[1,2-c]quinazoline derivatives: A novel class of potent and selective α_1 -adrenoceptor antagonists and antihypertensive agents. *J. Med. Chem.* **36**(15), 2196-2207 (1993).
2. Chern, J.-W., Shiau, C.-Y., and Lu, G.-Y. Studies on quinazolinones. 3: 1 novel and efficient route to the synthesis of conformationally restricted analogues of ketanserin and SGB-1534 as antihypertensive agents. *Bioorg. Med. Chem. Lett.* **1**(11), 571-574 (1991).

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, 04/30/2025

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