

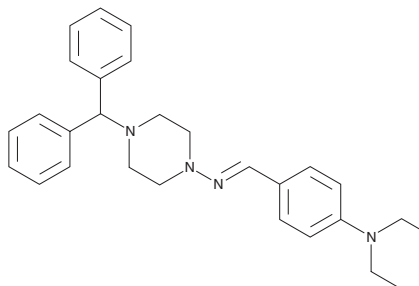
PRODUCT INFORMATION



L-138,037

Item No. 16526

CAS Registry No.: 31639-49-3
Formal Name: N-[[4-(diethylamino)phenyl]methylene]-4-(diphenylmethyl)-1-piperazinamine
MF: C₂₈H₃₄N₄
FW: 426.6
Purity: ≥98%
UV/Vis.: λ_{max}: 224, 318 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥2 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

L-138,037 is supplied as a crystalline solid. A stock solution may be made by dissolving the L-138,037 in the solvent of choice, which should be purged with an inert gas. L-138,037 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of L-138,037 in ethanol and DMSO is approximately 1 mg/ml and approximately 20 mg/ml in DMF.

Description

L-138,037 is a hydrazone derivative of 1-benzhydryl-4-aminopiperazine first studied for antimicrobial activity in the early 1970s. While this compound did not demonstrate *in vitro* antibacterial or antifungal activity, similar 1,4-disubstituted piperazines have been useful bactericides and fungicides demonstrating antimicrobial activity against *S. aureus*, *B. subtilis*, *C. perfringens*, *M. phlei*, and *S. cerevisiae* with MIC values ranging from 0.78-30 µg/ml.¹ L-138,037 has been explored for its ability to enhance binding of a ligand, such as FGF-2 or EPO, to a receptor, such as FGF-R1 or EPO-R, respectively, via non-covalent interaction with the ligand and with a second ligand (either a second molecule identical to the first ligand, or a second, different ligand molecule) for the same receptor, the receptor itself, or both.²

References

1. Yung, D.K., Mahony, D.E., and Whitehouse, L.W. Synthesis and *in vitro* antimicrobial evaluation of hydrazones of 1-phenyl-, 1-benzyl-, and 1-benzhydryl-4-aminopiperazines. *J. Pharm. Sci.* **60**(3), 386-389 (1971).
2. Soumya, R., Charette, M., and Finklestein, S.P. Methods and compositions for non-covalently enhanced receptor binding. *Biotrofix, Inc.* **WO2009/088975A2** (2009).

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, 09/08/2022

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