

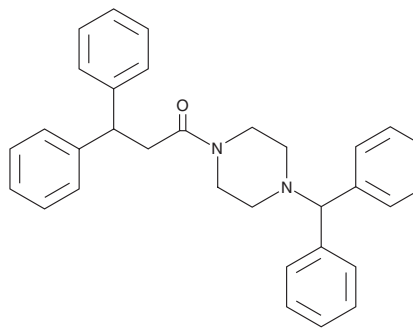
PRODUCT INFORMATION



NP-118809

Item No. 37539

CAS Registry No.: 41332-24-5
Formal Name: 1-[4-(diphenylmethyl)-1-piperazinyl]-3,3-diphenyl-1-propanone
Synonym: MK-6721
MF: C₃₂H₃₂N₂O
FW: 460.6
Purity: ≥95%
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

NP-118809 is supplied as a solid. A stock solution may be made by dissolving the NP-118809 in the solvent of choice, which should be purged with an inert gas. NP-118809 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of NP-118809 in these solvents is approximately 10 mg/ml.

Description

NP-118809 is an inhibitor of N-type voltage-gated calcium channels ($IC_{50} = 0.11 \mu M$).¹ It is selective for N-type over L-type voltage-gated calcium channels and human-ether-a-go-go (hERG), also known as K_v11.1 (IC_{50} s = 12.2 and 7.4 μM , respectively), as well as a panel of 112 other ion channels, receptors, and signaling targets at 10 μM . NP-118809 (25 mg/kg, i.p.) decreases the number of flinches in the second phase of the formalin test in a rat model of inflammatory pain.

Reference

1. Zamponi, G.W., Feng, Z.P., Zhang, L., *et al.* Scaffold-based design and synthesis of potent N-type calcium channel blockers. *Bioorg. Med. Chem. Lett.* **19**(22), 6467-6472 (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, 11/21/2023

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