

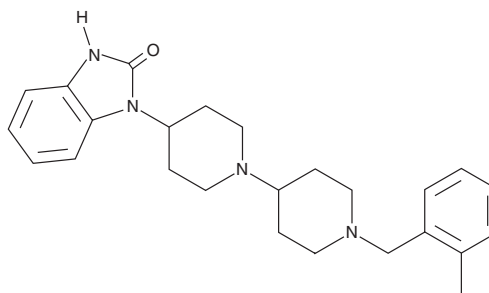
PRODUCT INFORMATION



TBPB

Item No. 22943

CAS Registry No.: 634616-95-8
Formal Name: 1,3-dihydro-1-[1'-[(2-methylphenyl)methyl][1,4'-bipiperidin]-4-yl]-2H-benzimidazol-2-one
MF: C₂₅H₃₂N₄O
FW: 404.6
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

TBPB is supplied as a crystalline solid. A stock solution may be made by dissolving the TBPB in the solvent of choice, which should be purged with an inert gas. TBPB is soluble in organic solvents such as chloroform and dimethyl formamide (DMF). The solubility of TBPB in these solvents is approximately 30 and 10 mg/ml, respectively.

TBPB is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, TBPB should first be dissolved in DMF and then diluted with the aqueous buffer of choice. TBPB has a solubility of approximately 0.25 mg/ml in a 1:3 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

TBPB is an allosteric agonist of M₁ muscarinic acetylcholine receptors (mAChRs; EC₅₀ = 158 nM in CHO-K1 cells expressing rat recombinant receptors).¹ It is selective for M₁ over M₂₋₅ receptors at concentrations up to 10 μM. TBPB (3 μM) potentiates NMDA-evoked currents in CA1 pyramidal cells. It increases production of the non-amyloidogenic amyloid precursor protein (APP) cleavage products APP_{sa} and CTFα in PC12 cells overexpressing the human M₁ receptor and APP when used at a concentration of 1 μM. TBPB decreases amphetamine-induced hyperlocomotion in a rat model that is predictive of antipsychotic-like activity.

References

1. Jones, C.K., Brady, A.E., Davis, A.A., et al. Novel selective allosteric activator of the M₁ muscarinic acetylcholine receptor regulates amyloid processing and produces antipsychotic-like activity in rats. *J. Neurosci.* **28(41)**, 10422-10433 (2008).

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, 11/28/2022

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