

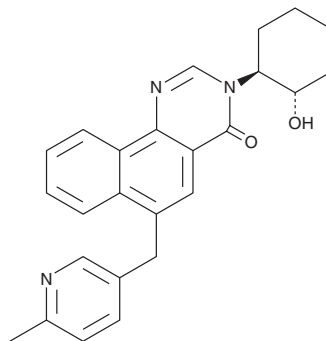
PRODUCT INFORMATION



MK-7622

Item No. 34407

CAS Registry No.: 1227923-29-6
Formal Name: 3-[(1S,2S)-2-hydroxycyclohexyl]-6-[(6-methyl-3-pyridinyl)methyl]-benzo[h]quinazolin-4(3H)-one
MF: C₂₅H₂₅N₃O₂
FW: 399.5
Purity: ≥98%
UV/Vis.: λ_{max}: 252 nm
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

MK-7622 is supplied as a solid. A stock solution may be made by dissolving the MK-7622 in the solvent of choice, which should be purged with an inert gas. MK-7622 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of MK-7622 in DMF is approximately 5 mg/ml and approximately 1 mg/ml in ethanol and DMSO.

MK-7622 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, MK-7622 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. MK-7622 has a solubility of approximately 0.2 mg/ml in a 1:4 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

MK-7622 is a positive allosteric modulator of muscarinic M₁ acetylcholine receptors (mAChRs).¹ It increases ACh-induced calcium flux in CHO cells expressing human M₁ receptors (EC₅₀ = 21 nM). MK-7622 (0.3 and 1 mg/kg) reverses scopolamine-induced cognitive deficits in an object retrieval detour task in rhesus macaques.

Reference

1. Uslaner, J.M., Kuduk, S.D., Wittmann, M., *et al.* Preclinical to human translational pharmacology of the novel M₁ positive allosteric modulator MK-7622. *J. Pharmacol. Exp. Ther.* **365**(3), 556-566 (2018).

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, 11/15/2022

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