

PRODUCT INFORMATION



Halopemide

Item No. 13205

CAS Registry No.: 59831-65-1
Formal Name: N-[2-[4-(5-chloro-2,3-dihydro-2-oxo-1H-benzimidazol-1-yl)-1-piperidinyl]-ethyl]-4-fluoro-benzamide

Synonyms: NSC 354856, R34301

MF: C₂₁H₂₂ClFN₄O₂

FW: 416.9

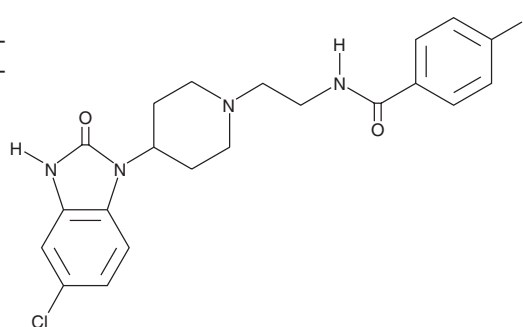
Purity: ≥98%

UV/Vis.: λ_{max}: 293 nm

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

Halopemide is supplied as a crystalline solid. A stock solution may be made by dissolving the halopemide in the solvent of choice, which should be purged with an inert gas. Halopemide is soluble in organic solvents such as DMSO and dimethyl formamide (DMF). The solubility of halopemide in these solvents is approximately 10 and 20 mg/ml, respectively.

Halopemide is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, halopemide should first be dissolved in DMF and then diluted with the aqueous buffer of choice. Halopemide has a solubility of approximately 0.1 mg/ml in a 1:9 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Halopemide is a potent inhibitor of phospholipase D (PLD), inhibiting human PLD₁ and PLD₂ *in vitro* (IC₅₀ = 220 and 310 nM, respectively) and PLD activity in cells.¹ Previously, halopemide has been found to inhibit dopamine receptors and was evaluated as a neuroleptic agent.²

References

1. Scott, S.A., Selvy, P.E., Buck, J.R., *et al.* Design of isoform-selective phospholipase D inhibitors that modulate cancer cell invasiveness. *Nat. Chem. Biol.* **5**(2), 108-117 (2009).
2. Seeman, P., Grigoriadis, D.E., and Niznik, H.B. Selectivity of agonists and antagonists at D₂ dopamine receptors compared to D₁ and S₂ receptors. *Drug Dev. Res.* **9**, 63-69 (1986).

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, 03/11/2024

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