

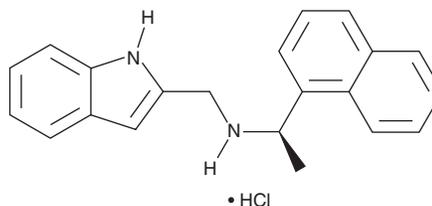
PRODUCT INFORMATION



Calindol (hydrochloride)

Item No. 17575

CAS Registry No.: 729610-18-8
Formal Name: N-[(1R)-1-(1-naphthalenyl)ethyl]-1H-indole-2-methanamine, monohydrochloride
MF: C₂₁H₂₀N₂ • HCl
FW: 336.9
Purity: ≥98%
UV/Vis.: λ_{max}: 224, 272, 282 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

Calindol (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the calindol (hydrochloride) in the solvent of choice, which should be purged with an inert gas. Calindol (hydrochloride) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of calindol (hydrochloride) in ethanol is approximately 15 mg/ml and approximately 50 mg/ml in DMSO and DMF.

Calindol (hydrochloride) is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, calindol (hydrochloride) should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Calindol (hydrochloride) has a solubility of approximately 0.3 mg/ml in a 1:2 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Calindol is a phenylalkylamine calcimimetic that facilitates the activation of the calcium-sensing receptor (CaSR), a G protein-coupled receptor that triggers phosphatidylinositol (PI)-intracellular calcium signaling and, in some cell types, cell proliferation.¹ It potently stimulates PI hydrolysis and proliferation through human CaSR (pEC₅₀ values are 6.9 and 7.4, respectively) in the presence of calcium.² Calindol directly interacts with CaSR in transmembrane domains and triggers PI hydrolysis less effectively in the absence of calcium, indicating that it is an allosteric modulator that cooperates with calcium to activate the CaSR.²⁻⁴ Calindol is used to study the properties and distribution of CaSR in different types of cells and tissues.^{4,5}

References

1. Kessler, A., Faure, H., Petrel, C., *et al. Bioorg. Med. Chem. Lett.* **14**(12), 3345-3349 (2004).
2. Ma, J.N., Owens, M., Gustafsson, M., *et al. J. Pharmacol. Exp. Ther.* **337**(1), 275-284 (2011).
3. Petrel, C., Kessler, A., Dauban, P., *et al. J. Biol. Chem.* **279**(18), 18990-18997 (2004).
4. Ray, K., Tisdale, J., Dodd, R.H., *et al. J. Biol. Chem.* **280**(44), 37013-37020 (2005).
5. Weston, A.H., Absi, M., Ward, D.T., *et al. Circ. Res.* **97**(4), 391-398 (2005).

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, 02/14/2024

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