

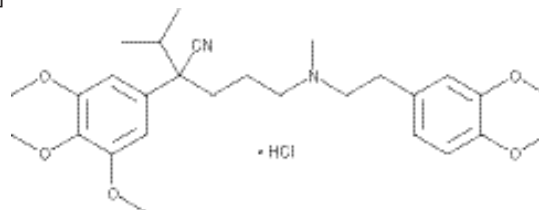
PRODUCT INFORMATION



(±)-Methoxyverapamil (hydrochloride)

Item No. 24270

CAS Registry No.: 16662-46-7
Formal Name: α-[3-[[2-(3,4-dimethoxyphenyl)ethyl]methylamino]propyl]-3,4,5-trimethoxy-α-(1-methylethyl)-benzeneacetonitrile, monohydrochloride
Synonyms: Gallopamil, NSC 274966
MF: C₂₈H₄₀N₂O₅ • HCl
FW: 521.1
Purity: ≥98%
UV/Vis.: λ_{max}: 281, 305 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

(±)-Methoxyverapamil (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the (±)-methoxyverapamil (hydrochloride) in the solvent of choice. (±)-Methoxyverapamil (hydrochloride) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of (±)-methoxyverapamil (hydrochloride) in ethanol and DMSO is approximately 10 mg/ml and approximately 16.7 mg/ml in DMF.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of (±)-methoxyverapamil (hydrochloride) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of (±)-methoxyverapamil (hydrochloride) in PBS, pH 7.2, is approximately 0.25 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

(±)-Methoxyverapamil is a phenylalkylamine L-type calcium channel blocker and methoxy derivative of verapamil (Item No. 14288).¹ (±)-Methoxyverapamil blocks recombinant rat L-type calcium channels expressed in tsA201 cells with IC₅₀ values of 782 and 8,000 nM at holding potentials of 10 and -60 mV, respectively. It blocks maitotoxin-induced calcium influx in NIH3T3 fibroblasts (IC₅₀ = 16 μM) and inhibits histamine-induced acid secretion in primary enriched guinea pig parietal cells (IC₅₀ = 10.9 μM).^{2,3} (±)-Methoxyverapamil binds to rat myocardial membranes (IC₅₀ = 16 nM) and inhibits electrically-stimulated contractions in rat right myocardial ventricular strips (EC₅₀ = 1.95 μM).⁴ In conscious, ovariectomized, post-partum rats, it decreases blood pressure by 30 mm Hg and induces 60% inhibition of uterine contractions when administered at doses of 204 and 324 μg/kg, respectively.⁵

References

1. Johnson, B.D., Hockerman, G.H., Scheuer, T., et al. *Mol. Pharmacol.* **50**(5), 1388-1400 (1996).
2. Daly, J.W., Lueders, J., Padgett, W.L., et al. *Biochem. Pharmacol.* **50**(8), 1187-1197 (1995).
3. Sewing, K.F. and Hannemann, H. *Pharmacology* **27**(1), 9-14 (1983).
4. Theodore, L.J., Nelson, W.L., Zobrist, R.H., et al. *J. Med. Chem.* **29**(9), 1789-1792 (1986).
5. Abel, M.H. and Hollingsworth, M. *Br. J. Pharmacol.* **85**(1), 263-269 (1985).

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, 03/19/2026

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