

PRODUCT INFORMATION



Mosapride (citrate)

Item No. 30048

CAS Registry No.: 112885-42-4

Formal Name: 4-amino-5-chloro-2-ethoxy-N-[[4-[(4-fluorophenyl)methyl]-2-morpholinyl]methyl]-benzamide, 2-hydroxy-1,2,3-monopropanetricarboxylate

Synonym: AS-4370

MF: $C_{21}H_{25}ClFN_3O_3 \cdot C_6H_8O_7$

FW: 614.0

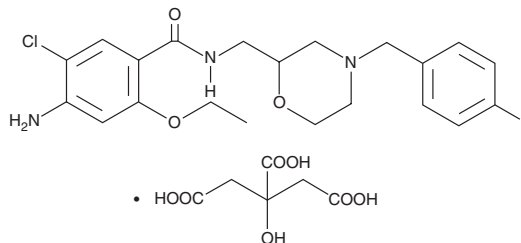
Purity: $\geq 98\%$

UV/Vis.: λ_{max} : 212, 232, 274, 309 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

Mosapride (citrate) is supplied as a solid. A stock solution may be made by dissolving the mosapride (citrate) in the solvent of choice, which should be purged with an inert gas. Mosapride (citrate) is soluble in organic solvents such as DMSO and dimethyl formamide (DMF). The solubility of mosapride (citrate) in these solvents is approximately 15 and 20 mg/ml, respectively.

Mosapride (citrate) is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, mosapride (citrate) should first be dissolved in DMF and then diluted with the aqueous buffer of choice. Mosapride (citrate) 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

Mosapride is an agonist of the serotonin (5-HT) receptor subtype 5-HT₄ ($K_i = 69.9$ nM).^{1,2} It is selective for 5-HT₄ over 5-HT₁, 5-HT₂, dopamine D₂, and α_1 - and α_2 -adrenergic receptors but acts as a partial antagonist at 5-HT₃ ($K_i = 1,189$ nM).^{1,2} It induces relaxation of precontracted rat esophageal thoracic muscularis mucosa preparations and enhances electrically evoked contractions in isolated guinea pig ileum (EC_{50} s = 208.4 and 73.2 nM, respectively).³ Mosapride (0.5 and 1 mg/kg, i.v.) increases antral and duodenal, but not ileal or colonic, motility in conscious dogs.¹ It also increases gastric emptying in rats when administered at doses ranging from 0.03 to 30 mg/kg.

References

1. Yoshida, N., Omoya, H., Oka, M., *et al.* AS-4370, a novel gastrokinetic agent free of dopamine D₂ receptor antagonist properties. *Arch. Int. Pharmacodyn. Ther.* **300**, 51-67 (1989).
2. Hirokawa, Y., Fujiwara, I., Suzuki, K., *et al.* Synthesis and structure-affinity relationships of novel N-(1-ethyl-4-methylhexahydro-1,4-diazepin-6-yl)pyridine-3-carboxamides with potent serotonin 5-HT₃ and dopamine D₂ receptor antagonistic activity. *J. Med. Chem.* **46**(5), 702-715 (2003).
3. Mine, Y., Yoshikawa, T., Oku, S., *et al.* Comparison of effect of mosapride citrate and existing 5-HT₄ receptor agonists on gastrointestinal motility *in vivo* and *in vitro*. *J. Pharmacol. Exp. Ther.* **283**(3), 1000-1008 (1997).

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, 01/03/2023

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