

PRODUCT INFORMATION

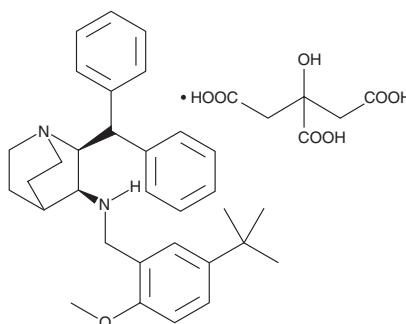


Maropitant (citrate)

Item No. 45625

CAS Registry No.: 862543-54-2
Formal Name: (2S,3S)-N-[[5-(1,1-dimethylethyl)-2-methoxyphenyl]methyl]-2-(diphenylmethyl)-1-azabicyclo[2.2.2]octan-3-amine, 2-hydroxy-1,2,3-propanetricarboxylate

MF: $C_{32}H_{40}N_2O \cdot C_6H_8O_7$
FW: 660.8
Purity: $\geq 98\%$
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

Maropitant (citrate) is supplied as a solid. A stock solution may be made by dissolving the maropitant (citrate) in the solvent of choice, which should be purged with an inert gas. Maropitant (citrate) is soluble (≥ 10 mg/ml) in ethanol and DMSO.

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 maropitant (citrate) can be prepared by directly dissolving the solid in aqueous buffers. Maropitant (citrate) is sparingly soluble (1-10 mg/ml) in PBS (pH 7.2). We do not recommend storing the aqueous solution for more than one day.

Description

Maropitant is a neurokinin-1 (NK_1) receptor antagonist.^{1,2} It reduces motion sickness- or doxorubicin-induced emesis in dogs. Formulations containing maropitant have been used as antiemetics in dogs.

References

1. Benchaoui, H.A., Siedek, E.M., De La Puente-Redondo, V.A., *et al.* Efficacy of maropitant for preventing vomiting associated with motion sickness in dogs. *Vet Rec.* **161**(13), 444-447 (2007).
2. Rau, S.E., Barber, L.G., and Burgess, K.E. Efficacy of maropitant in the prevention of delayed vomiting associated with administration of doxorubicin to dogs. *J. Vet. Intern. Med.* **24**(6), 1452-1457 (2021).

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, 07/31/2026

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