

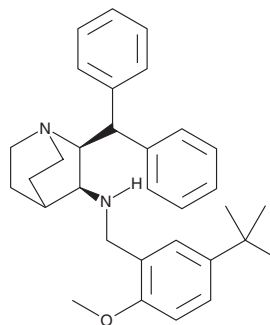
PRODUCT INFORMATION



Maropitant

Item No. 34816

CAS Registry No.: 147116-67-4
Formal Name: (2S,3S)-N-[[5-(1,1-dimethylethyl)-2-methoxyphenyl]methyl]-2-(diphenylmethyl)-1-azabicyclo[2.2.2]octan-3-amine
Synonym: CJ-11,972
MF: C₃₂H₄₀N₂O
FW: 468.7
Purity: ≥98%
UV/Vis.: λ_{max}: 277 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

Maropitant is supplied as a solid. A stock solution may be made by dissolving the maropitant in the solvent of choice, which should be purged with an inert gas. Maropitant is soluble in the organic solvent DMSO.

Description

Maropitant is an neurokinin-1 (NK₁) 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, 12/01/2022

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