

PRODUCT INFORMATION

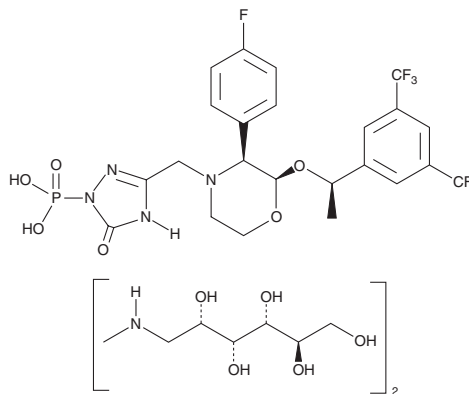


Fosaprepitant dimeglumine

Item No. 24034

CAS Registry No.: 265121-04-8
Formal Name: 1-deoxy-1-(methylamino)-D-glucitol, [3-[[[(2R,3S)-2-[[[(1R)-1-[3,5-bis(trifluoromethyl)phenyl]ethoxy]-3-(4-fluorophenyl)-4-morpholinyl]methyl]-2,5-dihydro-5-oxo-1H-1,2,4-triazol-1-yl]phosphonate (2:1)

Synonym: MK 0517
MF: C₂₃H₂₂F₇N₄O₆P • 2C₇H₁₇NO₅
FW: 1,004.8
Purity: ≥95%
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

Fosaprepitant dimeglumine is supplied as a crystalline solid. Aqueous solutions of fosaprepitant dimeglumine can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of fosaprepitant dimeglumine in PBS, pH 7.2, is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Fosaprepitant dimeglumine is a water-soluble prodrug form of the neurokinin-1 receptor antagonist aprepitant (Item No. 14867).¹ Formulations containing fosaprepitant dimeglumine have been used to prevent chemotherapy-induced nausea and vomiting.

Reference

1. Navari, R.M. Fosaprepitant (MK-0517): A neurokinin-1 receptor antagonist for the prevention of chemotherapy-induced nausea and vomiting. *Expert Opin. Investig. Drugs* **16(12)**, 1977-1985 (2007).

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, 11/10/2022

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