

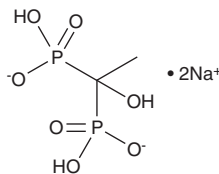
PRODUCT INFORMATION



Etidronate (sodium salt)

Item No. 20769

CAS Registry No.: 7414-83-7
Formal Name: P,P'-(1-hydroxyethylidene)bis-phosphonic acid, disodium salt
Synonyms: Etidronic Acid, Didronel, HEDP, Hydroxyethylidene Diphosphonic Acid
MF: C₂H₆O₇P₂ • 2Na
FW: 250.0
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

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

Description

Etidronate is a first generation, non-nitrogenous diphosphonate analog of endogenous pyrophosphate. It binds to hydroxyapatite and induces osteoclast apoptosis, thereby inhibiting abnormal bone resorption.¹ Unlike other bisphosphonates, etidronate may also prevent bone calcification.

Reference

1. Ohno, K., Mori, K., Orita, M., *et al.* Computational insights into binding of bisphosphates to farnesyl pyrophosphate synthase. *Curr. Med. Chem.* **18(2)**, 220-233 (2011).

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, 10/06/2022

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