

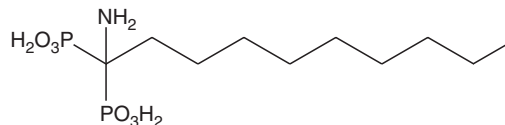
PRODUCT INFORMATION



1-Aminodecylidene *bis*-Phosphonic Acid

Item No. 13583

CAS Registry No.: 15049-93-1
Formal Name: (1-aminodecane-1,1-diyl)diphosphonic acid
Synonyms: 1-Aminodecane-1,1-Diphosphonic Acid, ARC 39
MF: C₁₀H₂₅NO₆P₂
FW: 317.3
Purity: ≥95%
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥2 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Description

1-Aminodecylidene *bis*-phosphonic acid is a potent inhibitor of acid sphingomyelinase (IC₅₀ = 20 nM).¹ It is selective for acid sphingomyelinase over neutral sphingomyelinase (IC₅₀ = >100 μM). 1-Aminodecylidene *bis*-phosphonic acid completely inhibits apoptosis induced by dexamethasone (Item No. 11015) in HepG2 cells.

Reference

1. Roth, A.G., Drescher, D., Yang, Y., *et al.* Potent and selective inhibition of acid sphingomyelinase by bisphosphonates. *Angew. Chem. Int. Ed. Engl.* **48(41)**, 7560-7563 (2009).

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, 02/09/2023

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