

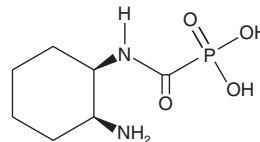
PRODUCT INFORMATION



cis-ACCP

Item No. 10012583

CAS Registry No.: 777075-44-2
Formal Name: P-[[[(1R,2S)-2-aminocyclohexyl]amino]carbonyl]-phosphonic acid
MF: C₇H₁₅N₂O₄P
FW: 222.2
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

cis-ACCP is supplied as a crystalline solid. A stock solution may be made by dissolving the cis-ACCP in water. The solubility of cis-ACCP in water is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Matrix metalloproteinases (MMPs) belong to a family of proteases that play a crucial role in tissue remodeling and repair by degrading extracellular matrix proteins, thus enabling cell migration.¹ Overexpression of MMPs are linked to diseases such as cancer, arthritis, osteoporosis, and arteriosclerosis. cis-ACCP is a reversible and competitive inhibitor of type IV collagen-specific MMP-2 and MMP-9 with preference towards MMP-2 (IC₅₀ = 4 and 20 μM, respectively).² At 100 μM, cis-ACCP prevents 90% of tumor cell invasion across matrigel-coated membranes *in vitro*.²

References

1. Nagase, H. and Woessner, J.F., Jr. Matrix metalloproteinases. *J. Biol. Chem.* **274**(31), 21491-21494 (1999).
2. Hoffman, A., Qadri, B., Frant, J., *et al.* Carbamoylphosphonate matrix metalloproteinase inhibitors 6: Cis-2-aminocyclohexylcarbamoylphosphonic acid, a novel orally active antimetastatic matrix metalloproteinase-2 selective inhibitor-synthesis and pharmacodynamic and pharmacokinetic analysis. *J. Med. Chem.* **51**, 1406-1414 (2008).

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, 04/08/2024

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