

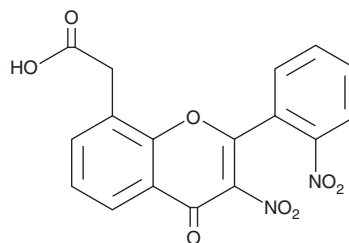
PRODUCT INFORMATION



Aminopeptidase N Inhibitor

Item No. 24307

CAS Registry No.: 596108-59-7
Formal Name: 3-nitro-2-(2-nitrophenyl)-4-oxo-4H-1-benzopyran-8-acetic acid
Synonym: AP-N Inhibitor
MF: C₁₇H₁₀N₂O₈
FW: 370.3
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.

Laboratory Procedures

Aminopeptidase N (AP-N) inhibitor is supplied as a crystalline solid. A stock solution may be made by dissolving the AP-N inhibitor in the solvent of choice. AP-N inhibitor is soluble in organic solvents such as DMSO and dimethyl formamide, which should be purged with an inert gas. The solubility of AP-N inhibitor in these solvents is approximately 33 mg/ml.

AP-N inhibitor is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, AP-N inhibitor should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. AP-N inhibitor has a solubility of approximately 0.2 mg/ml in a 1:4 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

AP-N inhibitor is a reversible inhibitor of AP-N/CD13 (IC₅₀ = 25 μM).¹ It is selective for AP-N/CD13 over matrix metalloproteinase-9 (MMP-9), angiotensin converting enzyme (ACE), neutral endopeptidase (NEP), γ-glutamyl transpeptidase, and the serine proteases dipeptidyl peptidase 4 (DPP-4) and cathepsin G at a concentration of 1 mM. AP-N inhibitor is non-cytotoxic to U937 cells at a concentration of 100 μM.

Reference

1. Bauvois, B., Puiffe, M.L., Bongui, J.B., *et al.* Synthesis and biological evaluation of novel flavone-8-acetic acid derivatives as reversible inhibitors of aminopeptidase N/CD13. *J. Med. Chem.* **46**(18), 3900-3913 (2003).

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, 03/28/2018

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