

PRODUCT INFORMATION



CP 471,474

Item No. 29442

CAS Registry No.: 210755-45-6
Formal Name: 2-[[[4-(4-fluorophenoxy)phenyl]sulfonyl]amino]-N-hydroxy-2-methyl-propanamide

MF: C₁₆H₁₇FN₂O₅S

FW: 368.4

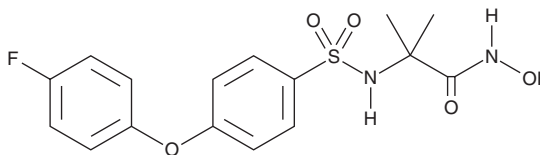
Purity: ≥98%

UV/Vis.: λ_{max}: 244 nm

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

CP 471,474 is supplied as a crystalline solid. A stock solution may be made by dissolving the CP 471,474 in the solvent of choice, which should be purged with an inert gas. CP 471,474 is soluble in the organic solvent DMSO.

Description

CP 471,474 is a broad-spectrum inhibitor of matrix metalloproteinases (MMPs) with IC₅₀ values of 0.7, 16, 13, and 0.9 nM for MMP-2, MMP-3, MMP-9, and MMP-13, respectively.¹ It is selective for these MMPs over MMP-1 (IC₅₀ = 1,170 nM). CP 471,474 reduces left ventricular enlargement and decreases the incidence of cardiac rupture in mouse models of myocardial infarction when administered at doses of 120 mg/kg twice per day and 240 mg/kg per day, respectively.^{1,2}

References

1. Rohde, L.E., Ducharme, A., Arroyo, L.H., *et al.* Matrix metalloproteinase inhibition attenuates early left ventricular enlargement after experimental myocardial infarction in mice. *Circulation* **99**(2), 3063-3070 (1999).
2. Fang, L., Gao, X.M., Moore, X.L., *et al.* Differences in inflammation, MMP activation and collagen damage account for gender difference in murine cardiac rupture following myocardial infarction. *J. Mol. Cell. Cardiol.* **43**(5), 535-544 (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, 11/11/2022

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