

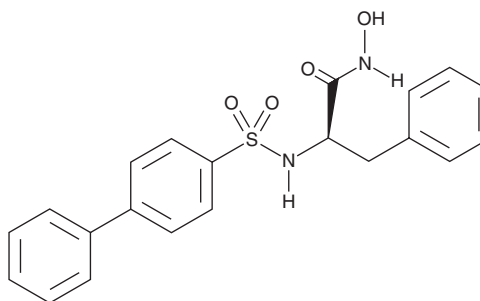
PRODUCT INFORMATION



MMP-2/MMP-9 Inhibitor II

Item No. 24314

CAS Registry No.: 193807-60-2
Formal Name: (αR)-α-[[[1,1'-biphenyl]-4-ylsulfonyl]amino]-N-hydroxy-benzenepropanamide
Synonym: Matrix Metalloproteinase-2/Matrix Metalloproteinase-9 Inhibitor II
MF: C₂₁H₂₀N₂O₄S
FW: 396.5
Purity: ≥98%
Supplied as: A 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

MMP-2/MMP-9 inhibitor II is supplied as a solid. A stock solution may be made by dissolving the MMP-2/MMP-9 inhibitor II in the solvent of choice. MMP-2/MMP-9 inhibitor II is soluble in organic solvents such as methanol and DMSO, which should be purged with an inert gas. The solubility of MMP-2/MMP-9 inhibitor II in these solvents is approximately 25 and 100 mg/ml, respectively.

Description

MMP-2/MMP-9 inhibitor II is a dual inhibitor of matrix metalloproteinase-2 (MMP-2) and MMP-9 (IC₅₀s = 17 and 30 nM, respectively).¹ In mice, it suppresses lung colonization of Lewis lung carcinoma cells and inhibits tumor-induced angiogenesis, tumor growth, and liver metastasis.^{1,2}

References

1. Tamura, Y., Watamane, F., Nakatani, T., *et al.* Highly selective and orally active inhibitors of type IV collagenase (MMP-9 and MMP-2): N-sulfonylamino acid derivatives. *J. Med. Chem.* **41**(4), 640-649 (1998).
2. Maki, H., Maekawa, R., Yoshida, H., *et al.* Antiangiogenic and antitumor effect of BPHA, an orally active matrix metalloproteinase inhibitor, in in vivo murine and human tumor model. *Gan To Kagaku Ryoho.* **26**(11), 1599-1606 (1999).

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, 10/05/2022

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