

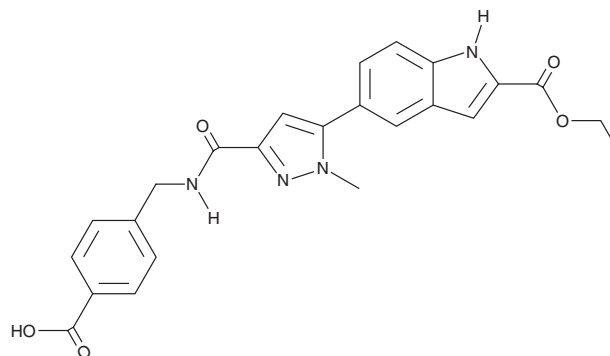
PRODUCT INFORMATION



BI-4394

Item No. 36855

CAS Registry No.: 1222173-37-6
Formal Name: 5-[3-[[[4-carboxyphenyl)methyl]amino]carbonyl]-1-methyl-1H-pyrazol-5-yl]-1H-indole-2-carboxylic acid, 2-ethyl ester
MF: C₂₄H₂₂N₄O₅
FW: 446.5
Purity: ≥95%
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

BI-4394 is supplied as a solid. A stock solution may be made by dissolving the BI-4394 in the solvent of choice, which should be purged with an inert gas. BI-4394 is soluble in the organic solvent DMSO at a concentration of approximately 10 mM.

Description

BI-4394 is an inhibitor of matrix metalloproteinase-13 (MMP-13; IC₅₀ = 0.001 μM for the recombinant human enzyme).¹ It is selective for MMP-13 over MMP-1-3, 7-10, -12, and -14 with IC₅₀ values ranging from 8.3 to greater than 22 μM for the recombinant human enzymes. BI-4394 inhibits the degradation of isolated bovine nasal cartilage explants cultured with human MMP-13 (IC₅₀ = 0.031 μM). It decreases disease severity in a mouse model of anti-type II collagen antibody and LPS-induced arthritis when administered at a dose of 100 mg/kg.²

References

1. Taylor, S.J., Abeywardane, A., Liang, S., *et al.* Fragment-based discovery of indole inhibitors of matrix metalloproteinase-13. *J. Med. Chem.* **54**(23), 8174-8187 (2011).
2. Taylor, S.J., Abeywardane, A., Liang, S., *et al.* Indole inhibitors of MMP-13 for arthritic disorders. *ACS Omega* **6**(29), 18635-18650 (2021).

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, 12/21/2022

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