

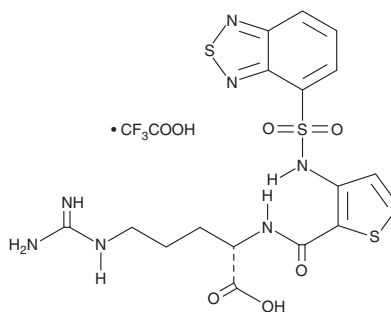
PRODUCT INFORMATION



EG 00229 (trifluoroacetate salt)

Item No. 33937

CAS Registry No.: 1210945-69-9
Formal Name: N²-[[3-[(2,1,3-benzothiazol-4-ylsulfonyl)amino]-2-thienyl]carbonyl]-L-arginine, 2,2,2-trifluoroacetate
MF: C₁₇H₁₉N₇O₅S₃ • CF₃COOH
FW: 611.6
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

EG 00229 (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the EG 00229 (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. EG 00229 (trifluoroacetate salt) is soluble in the organic solvent DMSO at a concentration of approximately 40 mg/ml.

Description

EG 00229 is an antagonist of neuropilin-1 (NRP-1).¹ It inhibits the binding of VEGF-A to porcine aortic endothelial cells expressing NRP-1, but not VEGFR2, with an IC₅₀ value of 8 μM. EG 00229 (100 μM) reduces VEGF-A-induced migration of human umbilical vein endothelial cells (HUVECs) and enhances the cytotoxicity of paclitaxel (Item No. 10461) or 5-fluorouracil (Item No. 14416) in A549 lung cancer cells. It also reduces severe acute respiratory syndrome coronavirus 2 (SARS-CoV-2) infectivity in Caco-2 cells when used at a concentration of 100 μM.²

References

1. Jarvis, A., Allerston, C.K., Jia, H., *et al.* Small molecule inhibitors of the neuropilin-1 vascular endothelial growth factor A (VEGF-A) interaction. *J. Med. Chem.* **53**(5), 2215-2226 (2010).
2. Daly, J.L., Simonetti, B., Klein, K., *et al.* Neuropilin-1 is a host factor for SARS-CoV-2 infection. *Science* **370**(6518), 861-865 (2020).

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, 10/31/2022

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