

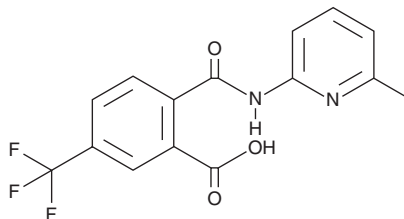
PRODUCT INFORMATION



AF38469

Item No. 35530

CAS Registry No.: 1531634-31-7
Formal Name: 2-[[[(6-methyl-2-pyridinyl)amino]carbonyl]-5-(trifluoromethyl)-benzoic acid
MF: C₁₅H₁₁F₃N₂O₃
FW: 324.3
Purity: ≥98%
UV/Vis.: λ_{max}: 283 nm
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

AF38469 is supplied as a solid. A stock solution may be made by dissolving the AF38469 in the solvent of choice, which should be purged with an inert gas. AF38469 is soluble in ethanol and DMSO.

Description

AF38469 is an inhibitor of sortilin, also known as the neurotensin-3 receptor (NTS₃; IC₅₀ = 330 nM).¹ It is selective for sortilin over neurotensin-1 (NTS₁), δ-opioid, angiotensin II type 1 (AT₁), endothelin A (ET_A), and ET_B receptors, GPR40, PPARδ, and matrix metalloproteinase-12 (MMP-12), and the prostaglandin E₂ (PGE₂) receptor subtype EP₁. AF38469 (400 nM) inhibits the epithelial-to-mesenchymal transition (EMT) in, and reduces migration and invasion of, U87 and A172 glioblastoma cells.² It also reduces tumor growth in a glioblastoma multiforme mouse xenograft model when administered at a dose of 0.01 g/kg per day.

References

1. Schröder, T.J., Christensen, S., Lindberg, S., *et al.* The identification of AF38469: An orally bioavailable inhibitor of the VPS10P family sorting receptor Sortilin. *Bioorg. Med. Chem. Lett.* **24**(1), 177-180 (2014).
2. Yang, W., Wu, P.-F., Ma, J.-X., *et al.* Sortilin promotes glioblastoma invasion and mesenchymal transition through GSK-3β/β-catenin/twist pathway. *Cell Death Dis.* **10**(3), 208 (2019).

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, 11/17/2022

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