

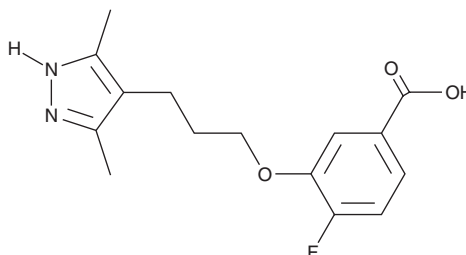
PRODUCT INFORMATION



Acoramidis

Item No. 41522

CAS Registry No.: 1446711-81-4
Formal Name: 3-[3-(3,5-dimethyl-1H-pyrazol-4-yl)propoxy]-4-fluoro-benzoic acid
Synonym: AG-10
MF: $C_{15}H_{17}FN_2O_3$
FW: 292.3
Purity: $\geq 98\%$
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

Acoramidis is supplied as a solid. A stock solution may be made by dissolving the acoramidis in the solvent of choice, which should be purged with an inert gas. Acoramidis is soluble (≥ 10 mg/ml) in DMSO and slightly soluble (0.1-1 mg/ml) in ethanol.

Description

Acoramidis is a kinetic stabilizer of transthyretin (TTR).¹ It selectively binds to wild-type TTR and TTR^{V122I} with negative cooperativity (K_{d1} s = 4.8 and 6.2 nM, respectively; K_{d2} s = 314 and 139 nM, respectively) over COX-1, COX-2, and the thyroid hormone nuclear receptor (IC_{50} s = >10 μM for all). Acoramidis inhibits pH-induced amyloidogenesis of wild-type TTR and TTR^{V122I}, indicating tetramer stabilization, when used at stoichiometric ratios of 1:1 and 1:0.5 with the number of TTR binding sites. It stabilizes TTR^{V122I} in serum isolated from patients with familial amyloid cardiomyopathy when used at a concentration of 10 μM .

Reference

1. Penchala, S.C., Connelly, S., Wang, Y., *et al.* AG10 inhibits amyloidogenesis and cellular toxicity of the familial amyloid cardiomyopathy-associated V122I transthyretin. *Proc. Natl. Acad. Sci. USA* **110**(24), 9992-9997 (2013).

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, 09/20/2024

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