

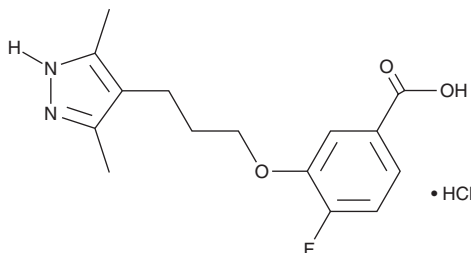
PRODUCT INFORMATION



Acoramidis (hydrochloride)

Item No. 41536

CAS Registry No.: 2242751-53-5
Formal Name: 3-[3-(3,5-dimethyl-1H-pyrazol-4-yl)propoxy]-4-fluoro-benzoic acid, monohydrochloride
Synonym: AG-10
MF: C₁₅H₁₇FN₂O₃ • HCl
FW: 328.8
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

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

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 μ M 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 μ M.

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, 09/16/2024

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