

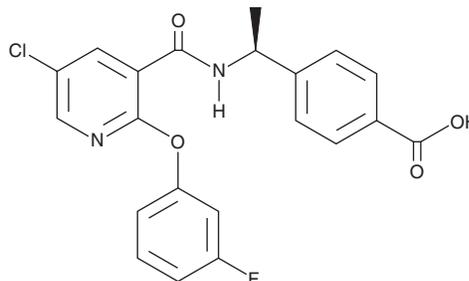
PRODUCT INFORMATION



AAT-008

Item No. 21810

CAS Registry No.: 847727-81-5
Formal Name: 4-[(1S)-1-[[[5-chloro-2-(3-fluorophenoxy)-3-pyridinyl]carbonyl]amino]ethyl]-benzoic acid
MF: C₂₁H₁₆ClFN₂O₄
FW: 414.8
Purity: ≥98%
UV/Vis.: λ_{max}: 234, 293 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥2 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

AAT-008 is supplied as a crystalline solid. A stock solution may be made by dissolving the AAT-008 in the solvent of choice. AAT-008 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of AAT-008 in ethanol is approximately 0.5 mg/ml and approximately 30 mg/ml in DMSO and DMF.

Description

AAT-008 is an orally bioavailable and potent antagonist of the prostaglandin E₂ (PGE₂) receptor subtype 4 (EP₄; IC₅₀ = 16.3 nM in a human EP₄ functional assay).¹ It is selective for EP₄ with IC₅₀ values of 2.4, 1,890, >20,000, and >20,000 nM for binding to human recombinant EP₄, EP₂, EP₁, and EP₃, respectively. AAT-008 has potent binding affinity for human, rat, and dog EP₄ (K_is = 0.97, 6.1, and 38 nM, respectively) and suppresses PGE₂-induced elevation of intracellular cAMP with an antagonistic potency (pA₂) of 1.1 nM *in vitro*. Oral administration of AAT-008 reduces carrageenan-induced mechanical hyperalgesia in rats in a dose-dependent manner.

Reference

1. Okumura, Y., Yamagishi, T., Nukui, S., *et al.* Discovery of AAT-008, a novel, potent, and selective prostaglandin EP₄ receptor antagonist. *Bioorg. Med. Chem. Lett.* **27(5)**, 1186-1192 (2017).

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, 07/31/2017

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