

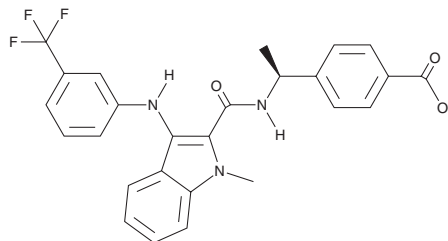
PRODUCT INFORMATION



AMX12006

Item No. 39027

CAS Registry No.: 2639775-01-0
Formal Name: 4-[(1S)-1-[[[1-methyl-3-[[3-(trifluoromethyl)phenyl]amino]-1H-indol-2-yl]carbonyl]amino]ethyl]-benzoic acid
MF: C₂₆H₂₂F₃N₃O₃
FW: 481.5
Purity: ≥98%
UV/Vis.: λ_{max}: 241 nm
Supplied as: A solid
Storage: -20°C
Stability: ≥3 years



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

Laboratory Procedures

AMX12006 is supplied as a solid. A stock solution may be made by dissolving the AMX12006 in the solvent of choice, which should be purged with an inert gas. AMX12006 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of AMX12006 in these solvents is approximately 1, 10, and 20 mg/ml, respectively.

Description

AMX12006 is an antagonist of the prostaglandin E₂ (PGE₂) receptor subtype EP₄ (IC₅₀ = 4.3 nM).¹ It is selective for EP₄ over EP₁, EP₂, and EP₃ receptors (IC₅₀s = >10 μM for all). AMX12006 is cytotoxic to MCF-7 and 4T1 breast and CT26 colon cancer cells (IC₅₀s = 46.73, 79.47, and 41.39 μM, respectively). It reduces tumor volume alone or in combination with the 5-fluorouracil prodrug capecitabine (Item No. 10487) in CT26 murine colon cancer models when administered at a dose of 150 mg/kg.

Reference

1. Das, D., Qiao, D., Liu, Z., *et al.* Discovery of novel, selective prostaglandin EP₄ receptor antagonists with efficacy in cancer models. *ACS Med. Chem. Lett.* **14**(6), 727-736 (2023).

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, 09/12/2023

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