

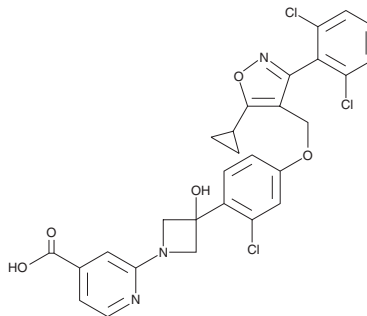
PRODUCT INFORMATION



Cilofexor

Item No. 25747

CAS Registry No.: 1418274-28-8
Formal Name: 2-[3-[2-chloro-4-[[5-cyclopropyl-3-(2,6-dichlorophenyl)-4-isoxazolyl]methoxy]phenyl]-3-hydroxy-1-azetidiny]-4-pyridinecarboxylic acid
MF: C₂₈H₂₂Cl₃N₃O₅
FW: 586.9
Purity: ≥98%
UV/Vis.: λ_{max}: 345 nm
Supplied as: A crystalline 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

Cilofexor is supplied as a crystalline solid. A stock solution may be made by dissolving the cilofexor in the solvent of choice, which should be purged with an inert gas. Cilofexor is soluble in organic solvents such as DMSO and dimethyl formamide (DMF). The solubility of cilofexor in these solvents is approximately 15 and 30 mg/ml, respectively.

Cilofexor is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, cilofexor should first be dissolved in DMF and then diluted with the aqueous buffer of choice. Cilofexor has a solubility of approximately 0.16 mg/ml in a 1:5 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Cilofexor inhibits binding of a synthetic peptide, comprising residues 676-700 of the steroid receptor coactivator 1 (SRC-1), to the farnesoid X receptor (FXR; EC₅₀ = <25 nM in a FRET activity assay).¹ It has FXR agonist activity in a mammalian one hybrid (M1H) assay (EC₅₀ = ≥100 nM).

Reference

1. Kinzel, O., Steeneck, C., and Kremoser, C. FXR (NR1H4) binding and activity modulating compounds. *Gilead Sciences Inc. US9539244B2* (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, 11/07/2022

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