

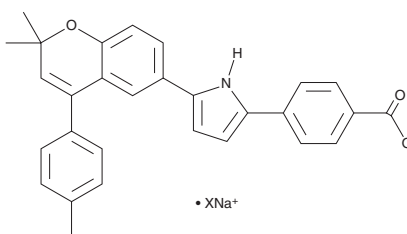
PRODUCT INFORMATION



YCT529

Item No. 39995

Formal Name: 4-(5-(2,2-dimethyl-4-(p-tolyl)-2H-chromen-6-yl)-1H-pyrrol-2-yl)benzoate, sodium salt
Synonym: GPHR-00354529
MF: C₂₉H₂₄NO₃ • XNa
FW: 434.5
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

YCT529 is supplied as a solid. A stock solution may be made by dissolving the YCT529 in the solvent of choice, which should be purged with an inert gas. YCT529 is soluble in DMSO. It is slightly soluble in acetonitrile.

YCT529 is slightly soluble in aqueous solutions. To enhance aqueous solubility, dilute the organic solvent solution into aqueous buffers or isotonic saline. If performing biological experiments, ensure the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

Description

YCT529 is an antagonist of retinoic acid receptor α (RAR α ; IC₅₀ = 6.8 nM).¹ It is selective for RAR α over RAR β and RAR γ (IC₅₀s = >3,700 nM for both). YCT529 (7.5 mg/kg per week) reversibly reduces epididymal sperm counts without reducing serum testosterone levels in male mice. It reduces germinal epithelial thickness in male dogs when administered at a dose of 25 mg/kg per day. YCT529 (10 mg/kg per week) decreases the number of embryos formed following copulation when administered to male rats.

Reference

1. Georg, G.I. and Cheryala, N. Contraceptive compounds and methods. *Regents of the University of Minnesota*, WO2022235693A1 (2022).

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, 03/07/2024

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