

PRODUCT INFORMATION



SKI2852

Item No. 45303

CAS Registry No.: 1346554-47-9

Formal Name: 2-[(2R)-4-[2-fluoro-4-(methylsulfonyl)phenyl]-2-methyl-1-piperazinyl]-N-[5-hydroxytricyclo[3.3.1.1.1]dec-2-yl]-4-pyrimidinecarboxamide

MF: C₂₇H₃₄FN₅O₄S

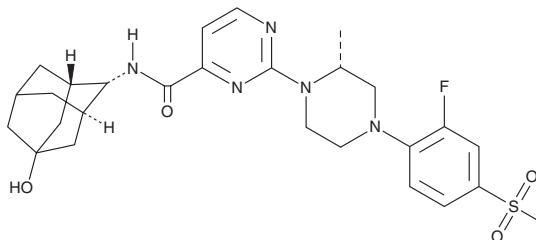
FW: 543.7

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

SKI2852 is supplied as a solid. A stock solution may be made by dissolving the SKI2852 in the solvent of choice, which should be purged with an inert gas. SKI2852 is soluble (≥10 mg/ml) in ethanol and DMSO.

Description

SKI2852 is an inhibitor of 11 β -hydroxysteroid dehydrogenase type 1 (11 β -HSD1; IC₅₀s = 2.9 and 1.6 nM for the human and mouse enzymes, respectively).¹ Oral administration of SKI2852 (20 mg/kg) reduces postprandial blood glucose, hemoglobin A1c (HbA1c), blood LDL and total cholesterol, and blood triglyceride levels in an *ob/ob* mouse model of type 2 diabetes and metabolic syndrome.

Reference

1. Ryu, J.H., Lee, J.A., Kim, S., *et al.* Discovery of 2-((R)-4-(2-fluoro-4-(methylsulfonyl)phenyl)-2-methylpiperazin-1-yl)-N-(((1R,2S,3S,5S,7S)-5-hydroxyadamantan-2-yl)pyrimidine-4-carboxamide (SKI2852): A highly potent, selective, and orally bioavailable inhibitor of 11 β -hydroxysteroid Dehydrogenase Type 1 (11 β -HSD1). *J. Med. Chem.* **59**(22), 10176-10189 (2016).

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, 03/09/2026

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