

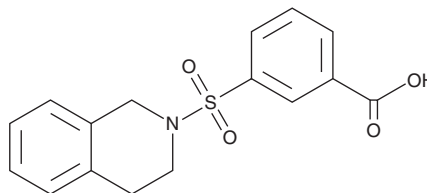
PRODUCT INFORMATION



CRT0036521

Item No. 36600

CAS Registry No.: 327092-81-9
Formal Name: 3-[(3,4-dihydro-2(1H)-isoquinolinyl)sulfonyl]-benzoic acid
MF: C₁₆H₁₅NO₄S
FW: 317.4
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

CRT0036521 is supplied as a solid. A stock solution may be made by dissolving the CRT0036521 in the solvent of choice, which should be purged with an inert gas. CRT0036521 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of CRT0036521 in these solvents is approximately 25 and 33 mg/ml, respectively. CRT0036521 is slightly soluble in ethanol.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of CRT0036521 can be prepared by directly dissolving the solid in aqueous buffers. The solubility of CRT0036521 in PBS (pH 7.2) is approximately 0.25 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

CRT0036521 is an inhibitor of prostaglandin F synthase, also known as aldo-keto reductase 1C3 (AKR1C3; IC₅₀ = 0.013 μM).¹ It is selective for prostaglandin F synthase/AKR1C3 over AKR1C1, AKR1C2, and AKR1C4 (IC₅₀s = 20, >30, and >30 μM, respectively).

Reference

1. Jamieson, S.M.F., Brooke, D.G., Heinrich, D., *et al.* 3-(3,4-Dihydroisoquinolin-2(1H)-ylsulfonyl)benzoic acids: Highly potent and selective inhibitors of the type 5 17-β-hydroxysteroid dehydrogenase AKR1C3. *J. Med. Chem.* **55**(17), 7746-7758 (2012).

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, 12/20/2022

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