

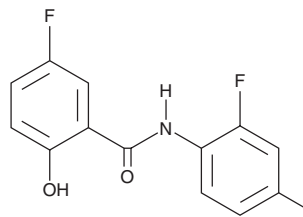
PRODUCT INFORMATION



NFATc1 Inhibitor A04

Item No. 41521

CAS Registry No.: 1912422-56-0
Formal Name: 5-fluoro-N-(2-fluoro-4-iodophenyl)-2-hydroxy-benzamide
Synonym: Nuclear Factor of Activated T cells c1 Inhibitor A04
MF: C₁₃H₈F₂INO₂
FW: 375.1
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

NFATc1 inhibitor A04 is supplied as a solid. A stock solution may be made by dissolving the NFATc1 inhibitor A04 in the solvent of choice, which should be purged with an inert gas. NFATc1 inhibitor A04 is soluble (≥10 mg/ml) in DMSO and sparingly soluble (1-10 mg/ml) in ethanol.

Description

NFATc1 inhibitor A04 is an inhibitor of nuclear factor of activated T cells c1 (NFATc1).¹ It inhibits RANKL-induced nuclear translocation of NFATc1 in RAW 264.7 cells when used at concentrations of 1.5 or 2 μM. NFATc1 inhibitor A04 reduces RANKL-induced differentiation of RAW 264.7 cells into osteoclasts (IC₅₀ = 1.57 μM).

Reference

1. Chen, C.L., Lee, C.C., Liu, F.L., *et al.* Design, synthesis and SARs of novel salicylanilides as potent inhibitors of RANKL-induced osteoclastogenesis and bone resorption. *Eur. J. Med. Chem.* **117**, 70-84 (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, 09/03/2024

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