

PRODUCT INFORMATION

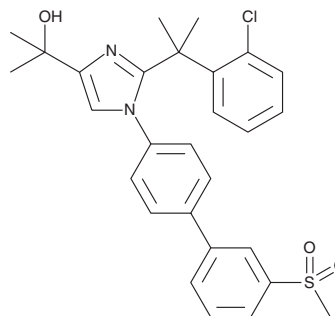


BMS 779788

Item No. 27339

CAS Registry No.: 918348-67-1
Formal Name: 2-[1-(2-chlorophenyl)-1-methylethyl]- α,α -dimethyl-1-[3'-(methylsulfonyl)[1,1'-biphenyl]-4-yl]-1H-imidazole-4-methanol

MF: C₂₈H₂₉ClN₂O₃S
FW: 509.1
Purity: \geq 98%
UV/Vis.: λ_{\max} : 258 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: \geq 4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

BMS 779788 is supplied as a crystalline solid. A stock solution may be made by dissolving the BMS 779788 in the solvent of choice, which should be purged with an inert gas. BMS 779788 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of BMS 779788 in these solvents is approximately 30 mg/ml.

Description

BMS 779788 is a partial agonist of liver X receptor α (LXR α) and LXR β ($K_{i,s}$ = 68 and 14 nM, respectively, in a radioligand binding assay).¹ It activates LXR α and LXR β with EC₅₀ values of 230 and 250 nM, respectively, in a transactivation assay. BMS 779788 increases expression of ATP-binding cassette transporter (ABCA1) in HeLa cells and isolated human and mouse whole blood (EC₅₀s = 33, 1,200, and 120 nM, respectively). It also increases expression of Abca1 and Abcg2 in isolated mouse blood cells but does not increase plasma or hepatic triglycerides when administered at a dose of 10 mg/kg.

References

1. Kick, E., Martin, R., Xie, Y., *et al.* Liver X receptor (LXR) partial agonists: Biaryl pyrazoles and imidazoles displaying a preference for LXR β . *Bioorg. Med. Chem. Lett.* **25**(2), 372-377 (2015).

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, 12/08/2022

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