

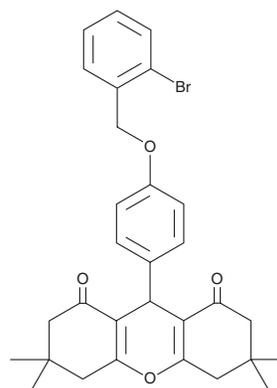
PRODUCT INFORMATION



BMS 986188

Item No. 21548

CAS Registry No.: 1776115-10-6
Formal Name: 9-[4-[(2-bromophenyl)methoxy]phenyl]-3,4,5,6,7,9-hexahydro-3,3,6,6-tetramethyl-1H-xanthene-1,8(2H)-dione
MF: C₃₀H₃₁BrO₄
FW: 535.5
Purity: ≥98%
UV/Vis.: λ_{max}: 223, 286 nm
Supplied as: A crystalline 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

BMS 986188 is supplied as a crystalline solid. A stock solution may be made by dissolving the BMS 986188 in the solvent of choice, which should be purged with an inert gas. BMS 986188 is soluble in the organic solvent dimethyl formamide (DMF). The solubility of BMS 986188 in DMF is approximately 1 mg/ml. BMS 986188 is also slightly soluble in ethanol and DMSO.

BMS 986188 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, BMS 986188 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. BMS 986188 has a solubility of approximately 0.3 mg/ml in a 1:2 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

BMS 986188 is a positive allosteric modulator of δ -opioid receptors, with an EC₅₀ value of 0.05 μ M in a β -arrestin recruitment assay in the presence, but not absence, of the opioid receptor agonist leu-enkephalin (Item No. 23283).¹ It is selective for δ - over μ -opioid receptors (EC₅₀ = >10 μ M) in the presence of leu-enkephalin and the μ -opioid receptor agonist endomorphin 1 (Item No. 23280), respectively.

Reference

1. Burford, N.T., Livingston, K.E., Canals, M., *et al.* Discovery, synthesis, and molecular pharmacology of selective positive allosteric modulators of the δ -opioid receptor. *J. Med Chem.* **58(10)**, 4220-4229 (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, 12/16/2022

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