

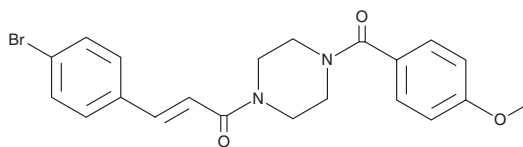
PRODUCT INFORMATION



NIBR189

Item No. 18792

CAS Registry No.: 1599432-08-2
Formal Name: 3-(4-bromophenyl)-1-[4-(4-methoxybenzoyl)-1-piperazinyl]-2E-propen-1-one
MF: C₂₁H₂₁BrN₂O₃
FW: 429.3
Purity: ≥95%
UV/Vis.: λ_{max}: 221, 282 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

NIBR189 is supplied as a crystalline solid. A stock solution may be made by dissolving the NIBR189 in the solvent of choice. NIBR189 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of NIBR189 in these solvents is approximately 1, 2, and 14 mg/ml, respectively.

NIBR189 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, NIBR189 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. NIBR189 has a solubility of approximately 0.1 mg/ml in a 1:8 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Epstein-Barr virus-induced G protein-coupled receptor 2 (EBI2, also known as GPR183) is an oxysterol receptor involved in B cell migration and function.¹ NIBR189 is an antagonist of EBI2 (IC₅₀s = 11 and 15 nM for human and mouse receptors, respectively).² It is selective for EBI2 over a panel of other receptors, enzymes, and transporters. NIBR189 blocks the oxysterol-induced migration of U937 monocytic cells.²

References

1. Sun, S. and Liu, C. 7α,25-dihydroxycholesterol-mediated activation of EBI2 in immune regulation and diseases. *Front. Pharmacol.* **6**, (2015).
2. Gessier, F., Preuss, I., Yin, H., et al. Identification and characterization of small molecule modulators of the Epstein-Barr virus-induced gene 2 (EBI2) receptor. *J. Med. Chem.* **57**, 3358-3368 (2014).

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, 10/07/2022

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