

# PRODUCT INFORMATION



## ML-401

Item No. 41450

**CAS Registry No.:** 1597489-14-9  
**Formal Name:** (2E)-3-(4-bromophenyl)-1-[4-

**MF:** C<sub>20</sub>H<sub>20</sub>BrClN<sub>2</sub>O

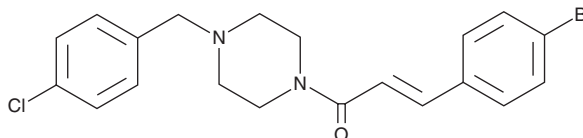
**FW:** 419.7

**Purity:** ≥98%

**Supplied as:** A solid

**Storage:** -20°C

**Stability:** ≥1 year



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

### Laboratory Procedures

ML-401 is supplied as a solid. A stock solution may be made by dissolving the ML-401 in the solvent of choice, which should be purged with an inert gas. ML-401 is slightly soluble (0.1-1 mg/ml) in DMSO and acetonitrile.

### Description

ML-401 is an Epstein-Barr virus-induced G protein-coupled receptor 2 (EBI2) antagonist (IC<sub>50</sub> = 1.03 nM).<sup>1</sup> It inhibits RS11846 cell chemotaxis (IC<sub>50</sub> = 4.533 nM).

### Reference

1. Ardecky, R., Sergienko, E., Zou, J., *et al.* Functional antagonists of EBI-2. (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, 10/22/2024

#### 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