

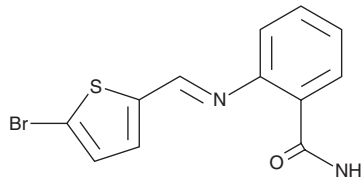
PRODUCT INFORMATION



BTYNB IMP1 Inhibitor

Item No. 25623

CAS Registry No.: 304456-62-0
Formal Name: 2-[[[(5-bromo-2-thienyl)methylene]amino]-benzamide
MF: C₁₂H₉BrN₂OS
FW: 309.2
Purity: ≥95%
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

BTYNB IMP1 inhibitor is supplied as a solid. A stock solution may be made by dissolving the BTYNB IMP1 inhibitor in the solvent of choice, which should be purged with an inert gas. BTYNB IMP1 inhibitor is soluble in the organic solvent DMSO.

Description

BTYNB IMP1 inhibitor is an inhibitor of the oncofetal mRNA-binding protein IMP1 (IC₅₀ = 5 μM for IMP1 binding to c-Myc mRNA).¹ It downregulates c-Myc and β-TrCP1 mRNA and reduces IMP1 protein levels in SK-MEL-2 and IGROV-1 cells. BTYNB IMP1 inhibits proliferation of IMP1-positive ES-2, IGROV-1, and SK-MEL-2 cells (IC₅₀s = 2.3, 3.6, and 4.5 μM, respectively) but not IMP1-negative BG-1 and T47D-KBluc cells (IC₅₀s = >50 μM for both). It also inhibits anchorage-independent growth of SK-MEL-2 cells when used at a concentration of 10 μM.

Reference

1. Mahapatra, L., Andruska, N., Mao, C., *et al.* A novel IMP1 inhibitor, BTYNB, targets c-Myc and inhibits melanoma and ovarian cancer cell proliferation. *Transl. Oncol.* **10(5)**, 818-827 (2017).

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, 12/07/2022

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