

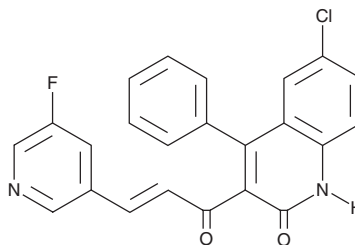
# PRODUCT INFORMATION



**SBI-0640756**

Item No. 40328

**Formal Name:** (E)-6-chloro-3-(3-(5-fluoropyridin-3-yl)acryloyl)-4-phenylquinolin-2(1H)-one  
**Synonym:** SBI-756  
**MF:** C<sub>23</sub>H<sub>14</sub>ClFN<sub>2</sub>O<sub>2</sub>  
**FW:** 404.8  
**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

SBI-0640756 is supplied as a solid. A stock solution may be made by dissolving the SBI-0640756 in the solvent of choice, which should be purged with an inert gas. SBI-0640756 is soluble in acetonitrile, methanol, and DMSO.

## Description

SBI-0640756 is an inhibitor of eukaryotic translation initiation factor 4 gamma 1 (eIF4G1).<sup>1</sup> It inhibits the association of eIF4G1 with the eIF4F complex in UACC-903 melanoma cell lysates in an m<sup>7</sup>GTP-agarose pull-down assay when used at a concentration of 0.75 μM. SBI-0640756 also inhibits eIF4G1-dependent and -independent mRNA translation in reporter assays using rabbit reticulocyte lysates (IC<sub>50</sub>s = 3.83 and 4.59 μM, respectively). It reduces the growth of 1205Lu, WM1346, A375, UACC-903, and WM3629 melanoma cells in a concentration-dependent manner. SBI-0640756 (0.5 mg/kg) decreases tumor incidence by 50% in an N-Ras<sup>Q61K</sup>-expressing, *Ink4a*<sup>-/-</sup> murine melanoma model. It also decreases tumor volume in an A375 mouse xenograft model when used in combination with the B-RAF<sup>V600E</sup> inhibitor PLX4720 (Item No. 15142).

## Reference

1. Feng, Y., Pinkerton, A.B., Hulea, L., *et al.* SBI-0640756 attenuates the growth of clinically unresponsive melanomas by disrupting the eIF4F translation initiation complex. *Cancer Res.* **75(24)**, 5211-5218 (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, 03/20/2024

## 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