

# PRODUCT INFORMATION



## Poloppin

Item No. 38330

CAS Registry No.: 683808-78-8

Formal Name: 5-(4-bromophenyl)-1-[2-(trifluoromethyl)phenyl]-1H-pyrrole-2-propanoic acid

Synonym: Polo Protein-protein Interaction Inhibitor

MF:  $C_{20}H_{15}BrF_3NO_2$

FW: 438.2

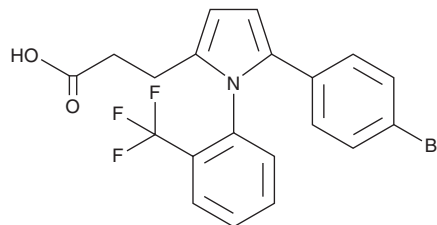
Purity:  $\geq 98\%$

UV/Vis.:  $\lambda_{max}$ : 295 nm

Supplied as: A solid

Storage:  $-20^\circ C$

Stability:  $\geq 4$  years



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

### Laboratory Procedures

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

### Description

Poloppin is an inhibitor of the protein-peptide interaction between polo-like kinase 1 (Plk1) and the Plk1-binding phosphopeptide ( $IC_{50} = 26.9 \mu M$ ).<sup>1</sup> It induces mitotic arrest in HeLa cells ( $EC_{50} = 29.9 \mu M$ ). Poloppin inhibits cell growth in murine pancreatic organoids expressing mutant K-Ras ( $GI_{50}s = 22.92-28.39 \mu M$ ).

### Reference

1. Narvaez, A.J., Ber, S., Crooks, A., *et al.* Modulating protein-protein interactions of the mitotic polo-like kinases to target mutant KRAS. *Cell Chem. Biol.* **24(8)**, 1017-1028 (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, 05/04/2023

#### 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