

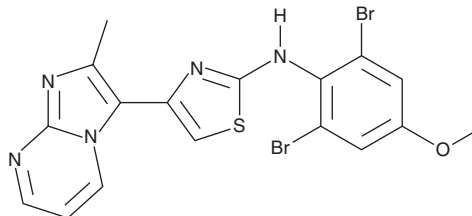
# PRODUCT INFORMATION



## PTC-209

Item No. 16277

**CAS Registry No.:** 315704-66-6  
**Formal Name:** N-(2,6-dibromo-4-methoxyphenyl)-4-(2-methylimidazo[1,2-a]pyrimidin-3-yl)-2-thiazolamine  
**MF:** C<sub>17</sub>H<sub>13</sub>Br<sub>2</sub>N<sub>5</sub>O<sub>2</sub>  
**FW:** 495.2  
**Purity:** ≥95%  
**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

PTC-209 is supplied as a crystalline solid. A stock solution may be made by dissolving the PTC-209 in the solvent of choice, which should be purged with an inert gas. PTC-209 is soluble in organic solvents such as DMSO and dimethyl formamide (DMF). The solubility of PTC-209 in these solvents is approximately 10 and 5 mg/ml, respectively.

PTC-209 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, PTC-209 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. PTC-209 has a solubility of approximately 0.1 mg/ml in a 1:7 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

Polycomb complex protein BMI-1 is a polycomb ring finger oncogene that regulates the p16 and p19 cell cycle inhibitor genes. It is necessary for efficient self-renewing cell divisions of stem cells in several tissues and can be over-expressed in tumors. PTC-209 is a BMI-1 inhibitor (IC<sub>50</sub> = ~ 0.5 μM) that irreversibly impairs colorectal cancer-initiating cell (CIC) growth.<sup>1</sup> It reduces tumor growth in CIC xenograft assays and abrogates colorectal cancer cell self-renewal *in vivo*, reducing their tumorigenic potential.<sup>1</sup>

### Reference

1. Kreso, A., van Galen, P., Pedley, N.M., *et al.* Self-renewal as a therapeutic target in human colorectal cancer. *Nat. Med.* **20**(1), 29-36 (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, 11/16/2022

#### 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