

# PTC-209

## BMI-1 inhibitor

PTC-209 is a potent and selective inhibitor of BMI-1 (B lymphoma Mo-MLV insertion region 1), with an IC50 value of 0.5  $\mu$ M. This compound is widely used in cancer research due to its ability to target cancer-initiating cells (CICs) and reduce tumor growth.

Key features and applications include:

- **Selective Inhibition:** Specifically targets BMI-1, a key regulator in the self-renewal of cancer stem cells.
- **Irreversible Action:** Impairs the growth and self-renewal properties of colorectal cancer-initiating cells.
- **High Potency:** Demonstrates significant anti-cancer activity in various cancer cell lines including colorectal, breast, lung, and cervical cancer cells.
- **Stem Cell Research:** Investigates the role of BMI-1 in the self-renewal and proliferation of cancer stem cells.
- **Combination Therapy:** Explored in combination with other anticancer drugs to enhance therapeutic efficacy.

Relevant disease states include:

- **Colorectal Cancer:** Reduces tumor growth and the self-renewal capacity of colorectal cancer-initiating cells.
- **Breast Cancer:** Inhibits the proliferation of breast cancer cells and reduces tumorigenic potential.
- **Lung Cancer:** Decreases the viability and growth of lung cancer cells.
- **Cervical Cancer:** Promotes cell cycle arrest and apoptosis in cervical cancer cell lines.

PTC-209 is a valuable tool in the field of oncology, providing insights into the mechanisms of cancer stem cell self-renewal and offering potential therapeutic strategies for various cancers.

- GMP format available

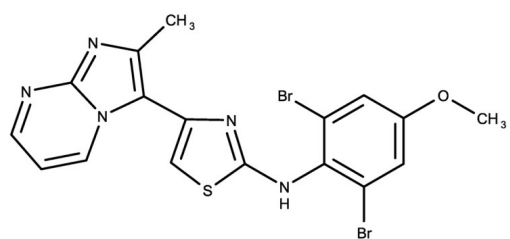
Ordering Information

Order Online »

ENZ-CHM354-0025	25mg
-----------------	------

Manuals, SDS & CofA

View Online »



## Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Solutions in DMSO may be stored at -20°C for up to 3 months.
Handling	Keep container tightly closed in a dry and well-ventilated place.
Short Term Storage	-20°C
Long Term Storage	-20°C
Shipping	Ambient Temperature

## Regulatory Status

RUO - Research Use Only

## Product Details

Alternative Name	N-(2,6-Dibromo-4-methoxyphenyl)-4-(2-methylimidazol[1,2-a]pyrimidin-3-yl)-2-thiazolamine
Appearance	Yellow solid.
CAS	315704-66-6
Couple Target	BMI-1
Couple Type	Inhibitor
Formula	$C_{17}H_{13}Br_2N_5OS$
Identity	Determined by NMR.
MW	495.19
Purity	≥98% HPLC
Solubility	Soluble in DMSO (>25 mg/ml).

Last modified: March 27, 2025



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-  
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-  
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-  
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-  
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)