

PTC-209 HBr

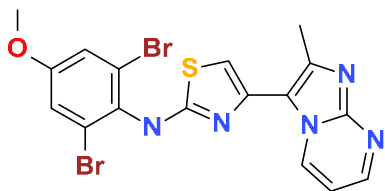
Cat. # SI9855

Background: Treatment of cells with PTC-209 leads to the selective reduction in the levels of BMI-1 protein (Kreso, van Galen et al. 2014). BMI-1 is a component of the polycomb PRC complex and interacts directly with the E3 ligase RING1A and/or RING2. The PRC complex mediates monoubiquitylation of histone H2A on Lys-119 and is involved in chromatin remodeling and repression of gene expression.

Application: Studying the function of BMI-1

Product Information

CAS No.:	1217022-63-3
Purity:	99%; NMR consistent with structure
Molecular Weight:	576.10
Physical State:	Powder
Quantity:	5mg
Solubility:	173mM in DMSO
Storage:	Store dessicated as supplied at -20°C for 2 years
Formula:	C ₁₇ H ₁₃ Br ₂ N ₅ OS.HBr
SMILES String:	Cc1c(n2cccnc2n1)c3csc(n3)Nc4c(cc(cc4Br)OC)Br

**References**

Kreso, A., P. van Galen, et al. (2014). "Self-renewal as a therapeutic target in human colorectal cancer." *Nat Med* **20**(1): 29-36.

All products are for research use only • not intended for human or animal diagnostic or therapeutic uses
Copyright © 2013 LifeSensors, Inc. All Rights Reserved