

## Product Information

<b>Description</b>	EPZ-5676 is an S-adenosyl methionine (SAM) competitive inhibitor of protein methyltransferase DOT1L with $K_i$ of 80 pM.
<b>Target:</b>	Histone Methyltransferase
<b>Pathway</b>	Chromatin/Epigenetics
<b>Chemical Name:</b>	(2R,3S,4S,5R)-2-(6-aminopurin-9-yl)-5-[[[3-[2-(6-tert-butyl-1H-benzimidazol-2-yl)ethyl]cyclobutyl]-propan-2-ylamino]methyl]oxolane-3,4-diol
<b>Molecular Formula:</b>	$C_{30}H_{42}N_8O_3$
<b>Supplied As:</b>	Solid
<b>Solubility:</b>	$\geq 28.15$ mg/ml in DMSO; insoluble in $H_2O$ ; $\geq 50.3$ mg/ml in EtOH with ultrasonic. For obtaining a higher solubility, please warm the tube at $37^\circ C$ and shake it in the ultrasonic bath for a while. Long-term storage in solution is not recommended. Not stable in diluted solutions.
<b>Stability:</b>	At least 6 months at $-20^\circ C$ . Avoid freeze/thaw cycles.
<b>Storage:</b>	$-20^\circ C$
<b>MW:</b>	562.71 Da
<b>Cas No.</b>	1380288-87-8
<b>Purity:</b>	$\geq 99\%$

## Chemical Structure

