

## Decitabine

Epigenetic Enzyme Inhibitor

### Catalog # D353-904B

Lot # P3320-15

CAS # 2353-33-5

### Product Description

Molecular Formula: C<sub>8</sub>H<sub>12</sub>N<sub>4</sub>O<sub>4</sub>

Appearance: White powder

Melting Point: 193-196°C

Molecular Weight: 228.2

Purity: >98%; NMR (Conforms)

Solubilization: May be dissolved in DMSO (25 mg/ml); or water (11 mg/ml). Ethanol (nil)

### Alias

5-Aza-2'-deoxycytidine; NSC-127716

### Specific Activity

DNA Methyl Transferase Inhibitor

### Storage and Stability

Store desiccated as supplied at 4°C for up to 2 years.

Store solutions at -20°C for up to 3 months.

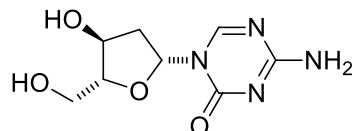
### Scientific Background

Inhibits DNA methyltransferase. Induces demethylation and reactivation of silenced genes suppressing the growth of human tumor cell lines<sup>1</sup>. Increases stem cell reprogramming efficiency. Stimulates osteogenic differentiation of human bone marrow stromal cells<sup>2</sup>. Induces differentiation and growth inhibition in IDH1 mutant glioma cells and decreases tumor growth in vivo<sup>3</sup>. Cell permeable.

### References

1. CM Bender et al. Cancer Res. 1998 58:95
2. AT El-Serafi et al. Differentiation 2011 81:35
3. S Turcan et al. Oncotarget 2013 4:1729

### Molecular Structure



## DECITABINE

Epigenetic Enzyme Inhibitor

Catalog # D353-904B

CAS # 2353-33-5

Lot # P3320-15

Purity >98%

Format White powder

Solubilization May be dissolved in DMSO (25 mg/ml); or water (11 mg/ml). Ethanol (nil)

Storage & Shipping Store desiccated as supplied at 4°C for up to 2 years. Store solutions at -20°C for up to 3 months.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: [orders@signalchem.com](mailto:orders@signalchem.com)  
[www.signalchem.com](http://www.signalchem.com)

**FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.**