

## 5-Azacytidine

Epigenetic Enzyme Inhibitor

**Catalog # D353-904**

Lot # P3320-2

CAS # 320-67-2

### Product Description

Molecular Formula: C<sub>8</sub>H<sub>12</sub>N<sub>4</sub>O<sub>5</sub>

Appearance: White powder

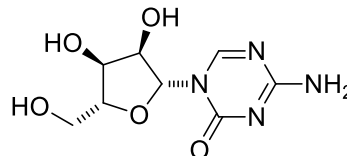
Melting Point: 226-232°C (dec.)

Molecular Weight: 244.2

Purity: >98% (TLC); NMR (Conforms)

Solubilization: May be dissolved in DMSO (25 mg/ml); or water (12 mg/ml).

### Molecular Structure



### Alias

4-Amino-1-(β-D-ribofuranosyl)-1,3,5-triazin-2(1H)-one,  
Ladakamycin

### Specific Activity

DNA hypomethylation agent

### Storage and Stability

Store desiccated as supplied at -20°C for up to 2 years.

Store solutions at -20°C for up to 1 month.

### Scientific Background

Specific inhibitor of DNA methyltransferase. Induces demethylation and reactivation of silenced genes. Increases stem cell reprogramming efficiency. Antitumor agent. Cell permeable.

### References

1. TS Mikkelsen et al. Nature 2008 454:49
2. AM Giraldo et al. Methods Mol. Biol. 2011 791:145

## 5-Azacytidine

Epigenetic Enzyme Inhibitor

Catalog # D353-904

CAS # 320-67-2

Lot # P3320-2

Purity >98%

Format White powder

Solubilization May be dissolved in DMSO (25 mg/ml); or water (12 mg/ml).

Storage & Shipping Store desiccated as supplied at -20°C for up to 2 years. Store solutions at -20°C for up to 1 month.

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.**