

---

## 5-Aza-2'-deoxycytidine

(Catalog No. M22000)

---

### Alternative Names

Decitabine; 5-Aza-2'-deoxy-D-cytidine

### Description

5-Aza-2'-deoxycytidine is an epigenetic modifier that inhibits DNA methyltransferase activity which results in DNA demethylation (hypomethylation) and gene activation by remodeling "opening" chromatin. Genes are synergistically reactivated when demethylation is combined with histone hyperacetylation.

### CAS Number

2353-33-5

### Chemical Formula

$C_8H_{12}N_4O_4$

### Molecular Weight

228.2

### Size

50 mg

### Purity

>95%

### Solubility

50% Acetic acid or DMSO

### Storage Temp

-20°C

### Ordering Information:

Products	Size	Cat. No.
5-Aza-2'-deoxycytidine	50 mg	M22000-1

*This product is for research purposes only. Not intended for use in diagnostic procedures.*