

---

## Zebularine

(Catalog No. M25001)

---

### Alternative Names

1-(β-D-Ribofuranosyl)-1,2-dihydropyrimidin-2-one, 2-Pyrimidone-1-β-D-ribose

### Description

A cell-permeable chemically stable cytidine analog that displays antitumor properties. Acts as transition state analog inhibitor of cytidine deaminase by binding at the active site as covalent hydrates. Also shown to inhibit DNA methylation and tumor growth both in vitro and in vivo.

### CAS Number

3690-10-6

### Chemical Formula

C<sub>9</sub>H<sub>12</sub>N<sub>2</sub>O<sub>5</sub>

### Molecular Weight

228.2

### Size

10 mg

### Purity

>95%

### Solubility

DMSO or MeOH

### Storage Temp

-20°C

### Ordering Information:

Products	Size	Cat. No.
Zebularine	10 mg	M25001 -1

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