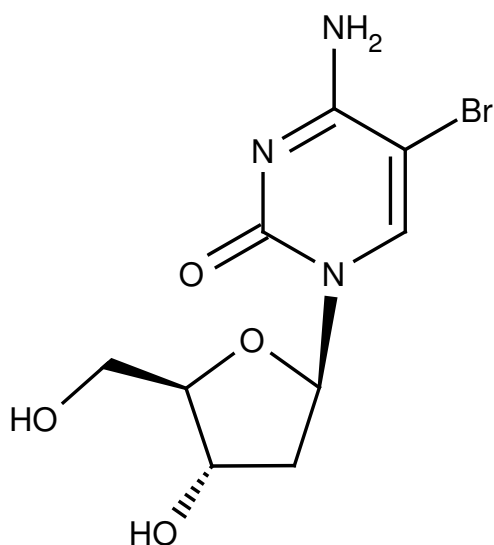


5-Bromo-2'-deoxy-cytidine

| Cat. No. | Amount |
|--------------|--------|
| N-DN-6496-1G | 1 g |
| N-DN-6496-5G | 5 g |



Molecular Formula: $C_9H_{12}BrN_3O_4$

Molecular Weight: 306.11 g/mol

Storage conditions: store at -20 °C

Shelf life: 30 months

Purity: $\geq 98\%$ (HPLC)

Appearance: white crystalline powder

Solubility: methanol

For research use only!