

5-Aza-2'-deoxycytidine (Decitabine)

Catalog No: 14100, 14101

Format: 50 mg, 10 mg

Chemical Properties:

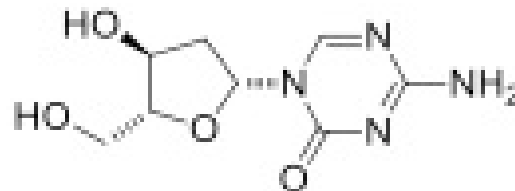
MW = 228.2

C₈H₁₂N₄O₄

CAS 2353-33-5

Physical Properties: White powder, MP = 193-196°C

Names: Decitabine, 5-Aza-2'-deoxycytidine, NSC-127716



Pharmacology: Inhibits DNA methyltransferases (DNMT). Induces demethylation and reactivation of silenced genes suppressing the growth of human tumor cell lines (ref 1). Increases stem cell reprogramming efficiency. Stimulates osteogenic differentiation of human bone marrow stromal cells (ref 2). Induces differentiation and growth inhibition in IDH1 mutant glioma cells and decreases tumor growth *in vivo* (ref 3). Cell permeable.

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

Fluorescent Properties: N/A

Quality Control:

>98% (TLC); NMR (Conforms)

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

Storage and Guarantee: Store desiccated as supplied at 4°C for up to 2 years. Store solutions at -20°C for up to 3 months. This product is guaranteed for 6 months from date of arrival.