

## 2'-Deoxy-5-fluorocytidine

**Catalog No:** 14108, 14109

**Format:** 250 mg, 50 mg

**Chemical Properties:**

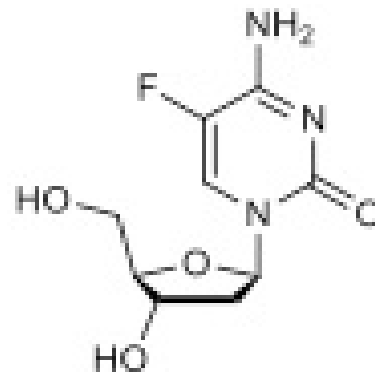
MW = 245.2

C<sub>9</sub>H<sub>12</sub>FN<sub>3</sub>O<sub>4</sub>

CAS 10356-76-0

Physical Properties: White solid; MP = 196°C

**Names:** 2'-Deoxy-5-fluorocytidine, 5-Fluoro-2'-deoxycytidine; FCDR; FdCyd; NSC-48006



**Pharmacology:** A DNA methyltransferase inhibitor currently in clinical trials for breast cancer and other solid tumors (ref 1). Like 5-Azacytidine and decitabine, 2'-Deoxy-5-fluorocytidine is a pyrimidine analog that integrates into chromatin to inhibit DNA methylation (ref 2). It blocks proliferation in colon cancer-derived HCT116 cells by activating DNA response pathways (ref 3). It inhibits various strains of pathogenic avian influenza viruses *in vitro* and *in vivo* (ref 4).

**Solubilization:** May be dissolved in DMSO (20 mg/ml)

**Fluorescent Properties:** N/A

**Quality Control:**

>98% (TLC); NMR (Conforms)

**References:**

1. J Ren *et al. Cell Signal.* 2011, 23:1082
2. H Gowher *et al. Cancer Biol. Ther.* 2004, 3:1062
3. Q Zhao *et al. Springerplus* 2002, 1:65
4. Y Kumaki *et al. Antiviral Res.* 2011, 92:329

**Storage and Guarantee:** Store desiccated as supplied at -20°C for up to 2 years. Store solutions at -20°C for up to 1 month. This product is guaranteed for 6 months from date of arrival.