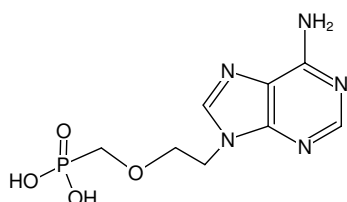


**Adefovir**

PMEA

P-[[2-(6-Amino-9H-purin-9-yl)ethoxy]methyl]phosphonic acid  
9-(2-phosphonylmethoxyethyl)adenine

Cat. No.	Amount
NU-281	100 mg



Structural formula of Adefovir

**For general laboratory use.****Shipping:** shipped at ambient temperature**Storage Conditions:** store at -20 °C

Short term exposure (up to 1 week cumulative) to ambient temperature possible.

**Shelf Life:** 12 months after date of delivery**Molecular Formula:** C<sub>8</sub>H<sub>12</sub>N<sub>5</sub>O<sub>4</sub>P (free acid)**Molecular Weight:** 273.19 g/mol (free acid)**Exact Mass:** 273.06 g/mol (free acid)**CAS#:** 106941-25-7**Purity:** ≥ 98 % (HPLC)**Form:** solid**Color:** white to off-white**Spectroscopic Properties:** λ<sub>max</sub> 259 nm, ε 15.4 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)