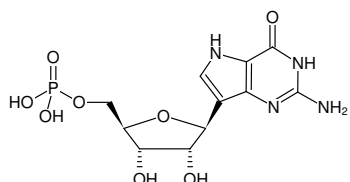


**9-Deaza-GMP**

9-Deazaguanosine-5'-monophosphate, Sodium salt

Cat. No.	Amount
NU-908S	100 µl (10 mM)
NU-908L	5 x 100 µl (10 mM)



Structural formula of 9-Deaza-GMP

**For research use only!****Shipping:** shipped on blue ice**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>11</sub>H<sub>15</sub>N<sub>4</sub>O<sub>8</sub>P (free acid)**Molecular Weight:** 362.23 g/mol (free acid)**Exact Mass:** 362.06 g/mol (free acid)**Purity:** ≥ 95 % (HPLC)**Form:** clear aqueous solution**Concentration:** 10 mM - 11 mM**pH:** 7.5 ±0.5**Spectroscopic Properties:** λ<sub>max</sub> 271 nm, ε 9.5 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)