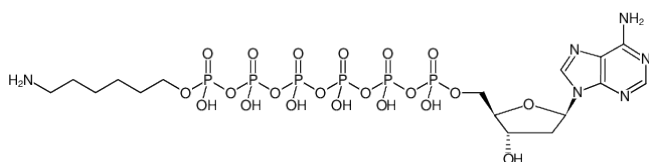



**zeta-(6-Aminoethyl)-dA6P**

zeta-(6-Aminoethyl)-2'-deoxyadenosine-5'-hexaphosphate, Sodium salt

Cat. No.	Amount
NU-282	50 µl (10 mM)



Structural formula of zeta-(6-Aminoethyl)-dA6P

**For general laboratory use.**
**Shipping:** shipped on gel packs

**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>16</sub>H<sub>32</sub>N<sub>6</sub>O<sub>21</sub>P<sub>6</sub> (free acid)

**Molecular Weight:** 830.3 g/mol (free acid)

**Exact Mass:** 830.0 g/mol (free acid)

**Purity:** ≥ 95 % (HPLC)

**Form:** solution in water

**Color:** colorless to slightly yellow

**Concentration:** 10 mM - 11 mM

**pH:** 7.5 ±0.5

**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)