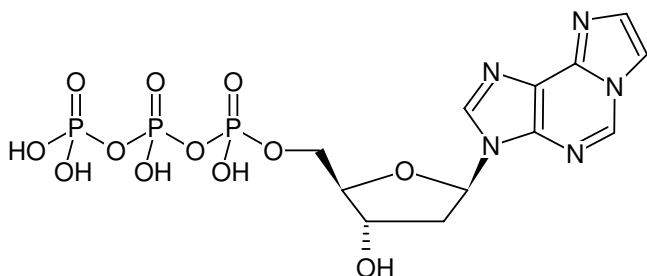


**Etheno-dATP (ϵ -dATP)**(1,N⁶-Etheno-dATP)1,N⁶-Etheno-2'-deoxyadenosine-5'-triphosphate, Sodium salt

Cat. No.	Amount
NU-256S	50 μ l (100 mM)
NU-256L	5 x 50 μ l (100 mM)

Selected References:

Revich *et al.* (1986) Utilization of 1,N⁶-Etheno-2'-deoxyadenosine-5'-triphosphate during DNA synthesis on natural templates, catalyzed by DNA polymerase I of *Escherichia coli*. *Carcinogenesis* **7 (9)**:1569.

Structural formula of Etheno-dATP (ϵ -dATP)**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:** C₁₂H₁₆N₅O₁₂P₃ (free acid)**Molecular Weight:** 515.20 g/mol (free acid)**Exact Mass:** 515.00 g/mol (free acid)**CAS#:** 81004-54-8**Purity:** \geq 95 % (HPLC)**Form:** solution in water**Color:** colorless to slightly yellow**Concentration:** 100 mM - 110 mM**pH:** 7.5 \pm 0.5**Spectroscopic Properties:** λ_{\max} 275 nm, ϵ 6.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5), λ_{exc} 300 nm, λ_{em} 415 nm