



Etheno-AppNHp (ϵ -AppNHp)

(1,N⁶-Etheno-AppNHp)

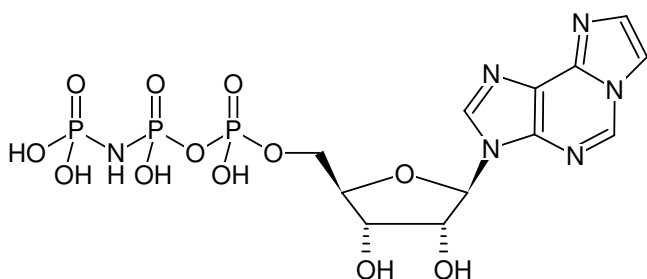
Etheno-AMPPNP (ϵ -AMPPNP)

1,N⁶-Etheno-adenosine-5'-[[β,γ]-imido]triphosphate, Sodium salt

Cat. No.	Amount
NU-1104	40 μ l (10 mM)

Selected References:

Yanagida (1981) Angles of nucleotides bound to cross-bridges in glycerinated muscle-fiber at various concentrations of Epsilon-ATP, Epsilon-ADP and Epsilon-AMPPNP detected by polarized fluorescence. *J. Mol. Biol.* **146** (4):539.



Structural formula of Etheno-AppNHp (ϵ -AppNHp)

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: 6 months after date of delivery

Molecular Formula: C₁₂H₁₇N₆O₁₂P₃ (free acid)

Molecular Weight: 530.22 g/mol (free acid)

Exact Mass: 530.01 g/mol (free acid)

CAS#: 78368-53-3

Purity: \geq 90 % (HPLC)

Form: solution in water

Color: colorless to slightly yellow

Concentration: 10 mM - 11 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