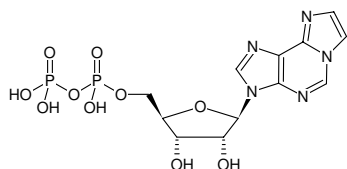



Etheno-ADP (ϵ -ADP)

 (1,N⁶-Etheno-ADP)

 1,N⁶-Etheno-adenosine-5'-diphosphate, Sodium salt

Cat. No.	Amount
NU-1140S	500 μ l (10 mM)
NU-1140L	5 x 500 μ l (10 mM)


 Structural formula of Etheno-ADP (ϵ -ADP)

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: C₁₂H₁₅N₅O₁₀P₂ (free acid)

Molecular Weight: 451.22 g/mol (free acid)

Exact Mass: 451.03 g/mol (free acid)

CAS#: 38806-39-2

Purity: \geq 95 % (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

Selected References:

 Anderson *et al.* (2011) A fluorescent GTP analog as a specific, high-precision label of microtubules. *Biotechniques*. **51** (1):43.

 Gräber *et al.* (1980) Phosphorylation of 1,N⁶-Etheno-ADP in Flash Groups and the concomitant Decay Kinetics of the Absorption Change at 515 nm. *Biochem. Biophys. Res. Comm.* **96** (3):1232.

 Secrist *et al.* (1972) Fluorescent modification of adenosine-containing coenzymes. Biological activities and spectroscopic properties. *Biochemistry* **11** (19):3499.