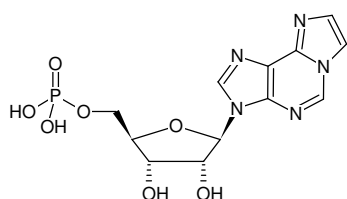


**Etheno-AMP (ϵ -AMP)**(1,N⁶-Etheno-AMP)1,N⁶-Etheno-adenosine-5'-monophosphate, Sodium salt

Cat. No.	Amount
NU-1141S	500 μ l (10 mM)
NU-1141L	5 x 500 μ l (10 mM)

Structural formula of Etheno-AMP (ϵ -AMP)**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:** C₁₂H₁₄N₅O₇P (free acid)**Molecular Weight:** 371.25 g/mol (free acid)**Exact Mass:** 371.06 g/mol (free acid)**CAS#:** 103213-41-8**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:**Latchezar *et al.* (2001) β_2 -Adrenoceptor-Mediated Prejunctional Facilitation and Postjunctional Inhibition of Sympathetic Neuroeffector Transmission in the Guinea Pig Vas Deferens. *J. Pharmacol. Exp. Ther.* **298**:623.Vandenbunner *et al.* (1976) 1,N⁶-Etheno-AMP and 1,N⁶-etheno-2'-deoxy-AMP as probes of the activator site of glycogen phosphorylase from rabbit skeletal muscle. *Proc. Nat. Acad. Sci. USA.* **73** (8):2696.