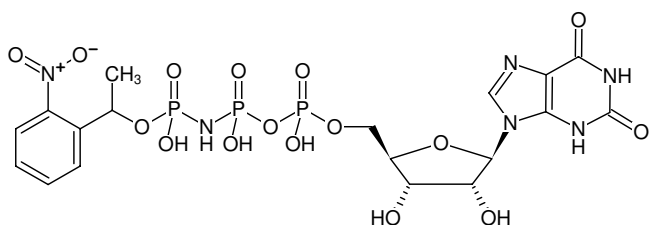


**NPE-caged-XppNHp**

(NPE-caged-XMPPNP)

Xanthosine-5'-[(β,γ)-imido]triphosphate, P³-(1-(2-nitrophenyl)-ethyl)-ester, Triethylammonium salt

Cat. No.	Amount
NU-308S	5 μ l (10 mM)
NU-308L	5 x 5 μ l (10 mM)



Structural formula of NPE-caged-XppNHp

For research use only!**Shipping:** shipped on gel packs**Storage Conditions:** store at -20 °C**Additional Storage Conditions:** store dark

Short term exposure (up to 1 week cumulative) to ambient temperature possible.

Shelf Life: 6 months after date of delivery**Molecular Formula:** C₁₈H₂₃N₆O₁₆P₃ (free acid)**Molecular Weight:** 672.33 g/mol (free acid)**Exact Mass:** 672.04 g/mol (free acid)**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} 254 nm, ϵ 13.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)