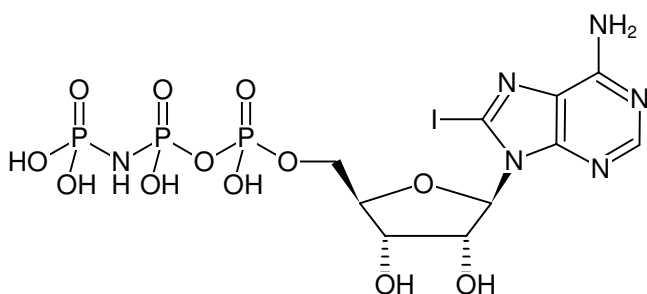


**8-Iodo-AppNHp**

(8I-AppNHp, 8I-AMPPNP)

8-Iodo-adenosine-5'-[(β,γ)-imido]triphosphate, Sodium salt

Cat. No.	Amount
NU-115	200 μ l (10 mM)



Structural formula of 8-Iodo-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:** C₁₀H₁₆N₆O₁₂P₃I (free acid)**Molecular Weight:** 632.09 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