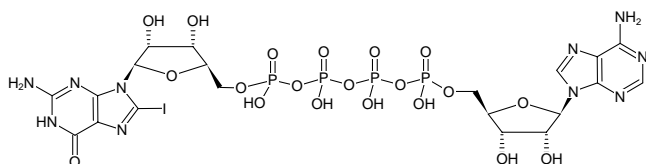


**AP₄(8-Iodo-G)**(AP₄(8I-G))P¹-(5'-Adenosyl) P⁴-[5'-(8-iodo)-guanosyl] tetraphosphate, Sodium salt

Cat. No.	Amount
NU-510S	30 µl (10 mM)
NU-510L	5 x 30 µl (10 mM)

Structural formula of AP₄(8-Iodo-G)**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:** C₂₀H₂₇IN₁₀O₂₀P₄ (free acid)**Molecular Weight:** 978.28 g/mol (free acid)**Purity:** ≥ 95 % (HPLC)**Form:** clear aqueous solution**Concentration:** 10 mM - 11 mM**pH:** 7.5 ±0.5**Spectroscopic Properties:** λ_{max} 259 nm; ε 27.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)