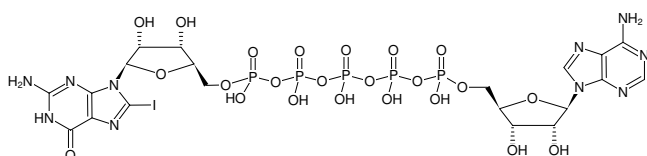


**AP₅(8-Iodo-G)**(AP₅(8I-G))P¹-(5'-Adenosyl) P⁵-[5'-(8-iodo)-guanosyl] pentaphosphate, Sodium salt

| Cat. No. | Amount |
|----------|-------------------|
| NU-511S | 30 µl (10 mM) |
| NU-511L | 5 x 30 µl (10 mM) |

Structural formula of AP₅(8-Iodo-G)**For research use only!****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:** C₂₀H₂₈IN₁₀O₂₃P₅ (free acid)**Molecular Weight:** 1058.26 g/mol (free acid)**Exact Mass:** 1057.91 g/mol (free acid)**Purity:** ≥ 95 % (HPLC)**Form:** solution in water**Color:** colorless to slightly yellow**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)