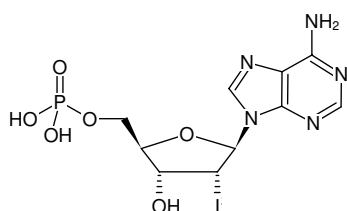


**2'-Iodo-dAMP**

(2'-I-dAMP)

2'-Iodo-2'-deoxyadenosine-5'-monophosphate, Sodium salt

Cat. No.	Amount
NU-147	250 µl (10 mM)



Structural formula of 2'-Iodo-dAMP

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:** C₁₀H₁₃N₅O₆PI (free acid)**Molecular Weight:** 457.12 g/mol (free acid)**Exact Mass:** 456.96 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 15.1 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.0)