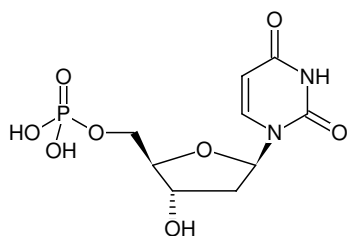


**dUMP**

2'-Deoxyuridine-5'-monophosphate, Disodium salt

Cat. No.	Amount
NU-1030S	500 µl (10 mM)
NU-1030L	5 x 500 µl (10 mM)



Structural formula of dUMP

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:** C₉H₁₃N₂O₈P (free acid)**Molecular Weight:** 308.18 g/mol (free acid)**CAS#:** 42155-08-8**EC number:** 255-687-1**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} 262 nm, ε 10.0 L mmol⁻¹ cm⁻¹
(Tris-HCl pH 7.5)