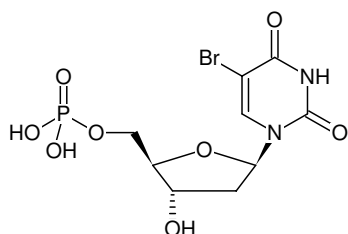


**5-Bromo-dUMP**

(5Br-dUMP)

5-Bromo-2'-deoxyuridine-5'-monophosphate, Sodium salt

Cat. No.	Amount
NU-1031S	100 µl (10 mM)
NU-1031L	5 x 100 µl (10 mM)



Structural formula of 5-Bromo-dUMP

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: 12 months after date of delivery**Molecular Formula:** C₉H₁₂N₂O₈PBr (free acid)**Molecular Weight:** 387.08 g/mol (free acid)**Exact Mass:** 385.95 g/mol (free acid)**CAS#:** 51432-32-7**EC number:** 257-200-8**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} 278 nm, ε 9.7 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)