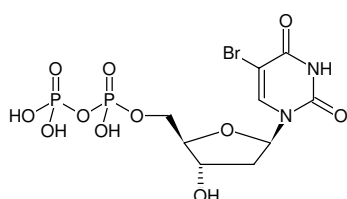


**5-Bromo-dUDP**

(5Br-dUDP)

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

Cat. No.	Amount
NU-129S	30 µl (10 mM)
NU-129L	5 x 30 µl (10 mM)



Structural formula of 5-Bromo-dUDP

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:** C₉H₁₃N₂O₁₁P₂Br (free acid)**Molecular Weight:** 467.06 g/mol (free acid)**Exact Mass:** 465.92 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} 278 nm, ε 9.7 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)