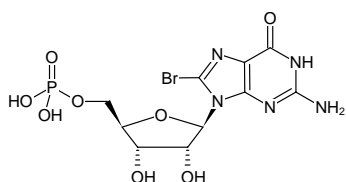


**8-Bromo-GMP**

(8Br-GMP)

8-Bromo-guanosine-5'-monophosphate, Sodium salt

Cat. No.	Amount
NU-144S	100 µl (10 mM)
NU-144L	5 x 100 µl (10 mM)



Structural formula of 8-Bromo-GMP

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:** C₁₀H₁₃N₅O₈PBr (free acid)**Molecular Weight:** 442.11 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 14.0 L mmol⁻¹ cm⁻¹
(Tris-HCl pH 7.5)