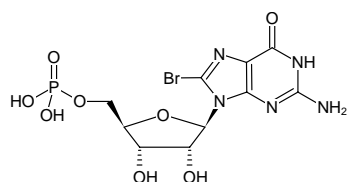


**8-Bromo-GMP**

(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. If stored as recommended, Jena Bioscience guarantees optimal performance of this product for 12 months after date of delivery.

**Shelf Life:** 12 months**Molecular Formula:** C<sub>10</sub>H<sub>13</sub>N<sub>5</sub>O<sub>8</sub>PBr (free acid)**Molecular Weight:** 442.11 g/mol (free acid)**Purity:** ≥ 95 % (HPLC)**Form:** clear aqueous solution, pH 7.5 ±0.5**Concentration:** 10 mM - 11 mM**pH:** 7.5 ±0.5

**Spectroscopic Properties:** λ<sub>max</sub> 259 nm; ε 14.0 L mmol<sup>-1</sup> cm<sup>-1</sup>  
(Tris-HCl pH 7.5)