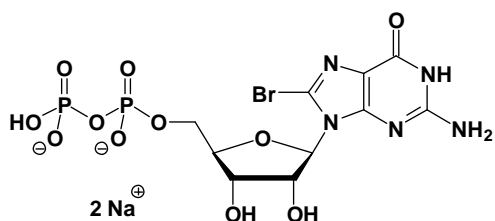


8Br-GDP

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

Cat. No.	Amount
NU-124S	50 Units
NU-124L	250 Units



Cat. No.: NU-124

Molecular Formula: $C_{10}H_{12}N_5O_{11}P_2Br$ (Anion)

Molecular Weight: 520.08 (Anion)

Purity: > 95%, clear aqueous solution, pH 7.5

Storage conditions:

Short term exposure (up to 1 week cumulative) to ambient temperature possible. Long term storage at $< -20^{\circ}C$. If stored as recommended, Jena Bioscience guarantees optimal performance of this product for 12 months after date of delivery.

For research use only!

1 unit = 1 μ l of a 10 mM solution