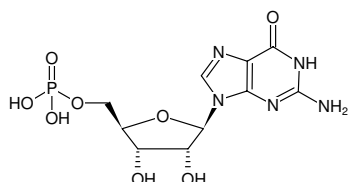


**GMP - Solid**

Guanosine-5'-monophosphate, Sodium salt

Cat. No.	Amount
NU-1028-5G	5 g
NU-1028-25G	25 g



Structural formula of GMP - Solid

**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:** C<sub>10</sub>H<sub>14</sub>N<sub>5</sub>O<sub>8</sub>P (free acid)**Molecular Weight:** 363.22 g/mol (free acid)**Exact Mass:** 363.06 g/mol (free acid)**CAS#:** 5550-12-9**Purity:** ≥ 95 % (HPLC)**Form:** solid**Color:** white to off-white**Spectroscopic Properties:** λ<sub>max</sub> 252 nm, ε 13.7 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)