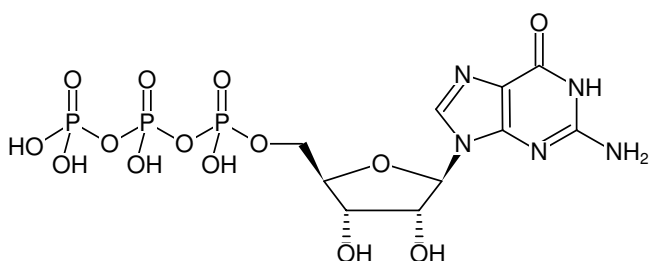


**GTP - Solid**

Guanosine 5'-triphosphate, Sodium salt

Cat. No.	Amount
NU-1047-200	200 mg
NU-1047-1G	1 g
NU-1047-10G	10 g
NU-1047-100G	100 g

**Quality Control Specifications:** in vitro transcription (T7 RNA polymerase): visible RNA fragments after 5 min incubation, Proteases: not detectable



Structural formula of GTP - Solid

**For general laboratory use.**

**Shipping:** Shipped on gel packs

**Storage Conditions:** Store at -20 °C

**Shelf Life:** 12 months

**Molecular Formula:** C<sub>10</sub>H<sub>16</sub>N<sub>5</sub>O<sub>14</sub>P<sub>3</sub> (free acid)

**Molecular Weight:** 523.18 g/mol (free acid)

**Exact Mass:** 522.99 g/mol (free acid)

**CAS#:** 36051-31-7

**Purity:** ≥ 90 % (HPLC)

**Form:** Powder

**Color:** White to slightly yellow

**Solubility:** Soluble in water. Clear solution.

**Spectroscopic Properties:** λ<sub>max</sub> 252 nm, ε 14.2 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.0)