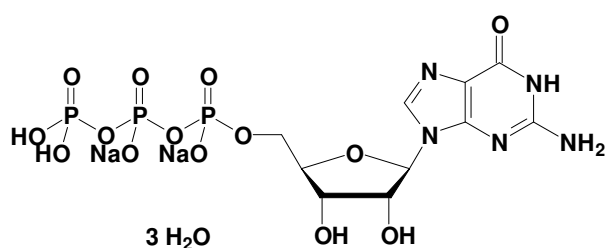


GTP, lyophilized

Guanosine 5'-triphosphate, sodium salt

Cat. No.	Amount
NU-1012-100	100 mg, lyophilized
NU-1012-1G	1 g, lyophilized
NU-1012-10G	10 g, lyophilized
NU-1012-100G	100 g, lyophilized



Molecular Formula: $C_{10}H_{14}N_5Na_2O_{14}P_3 \times 3H_2O$

Molecular Weight: 621.18

Purity: >90%

Storage conditions:

Short term exposure (up to 1 week cumulative) to ambient temperature possible. Long term storage at < -20°C.

For research use only!