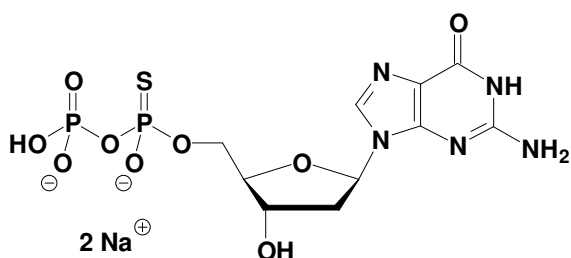


## dGDP $\alpha$ S

2'-Deoxy-Guanosine-5'-( $\alpha$ -thio)-diphosphate, Sodium salt

Cat. No.	Amount
NU-1164S	20 Units
NU-1164L	100 Units



**Cat. No.:** NU-1164

**Molecular Formula:** C<sub>10</sub>H<sub>13</sub>N<sub>5</sub>O<sub>9</sub>P<sub>2</sub>S (Anion)

**Molecular Weight:** 441.24 (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°C. If stored as recommended, Jena Bioscience guarantees optimal performance of this product for 12 months after date of delivery.

\* 1 unit = 1  $\mu$ l of a 10 mM solution

**For research use only!**