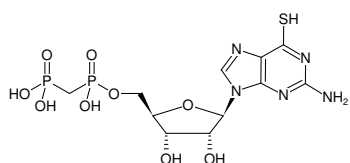


**6-Thio-GpCp**

(6-Thio-GMPCP)

6-Thio-guanosine-5'-[( $\alpha,\beta$ )-methylene]diphosphate, Sodium salt

Cat. No.	Amount
NU-417S	30 $\mu$ l (10 mM)
NU-417L	5 x 30 $\mu$ l (10 mM)



Structural formula of 6-Thio-GpCp

**For research use only!****Shipping:** shipped on gel packs**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>11</sub>H<sub>17</sub>N<sub>5</sub>O<sub>9</sub>P<sub>2</sub>S (free acid)**Molecular Weight:** 457.29 g/mol (free acid)**Exact Mass:** 457.02 g/mol (free acid)**Purity:**  $\geq$  95 % (HPLC)**Form:** solution in water**Color:** colorless to slightly yellow**Concentration:** 10 mM - 11 mM**pH:** 7.5  $\pm$  0.5**Spectroscopic Properties:**  $\lambda_{\max}$  257 nm,  $\epsilon$  8.8 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5),  $\lambda_{\max}$  342 nm,  $\epsilon$  24.8 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)