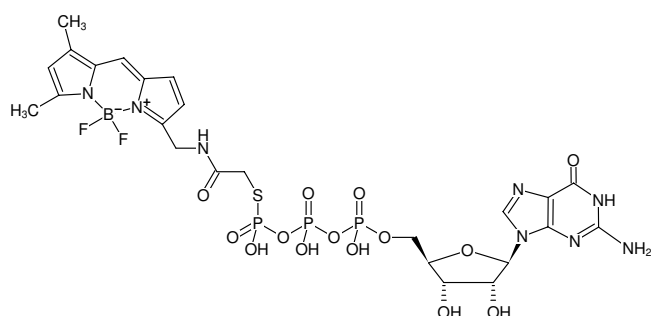


**GTPyS-BDP-FL**

also known as BODIPY® FL GTPyS

Guanosine-5'-(γ-thio)-triphosphate, labeled with BODIPY® FL, Sodium salt

Cat. No.	Amount
NU-973	50 µl (5 mM)



Structural formula of GTPyS-BDP-FL

**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>24</sub>H<sub>30</sub>BF<sub>2</sub>N<sub>8</sub>O<sub>14</sub>P<sub>3</sub>S (free acid)**Molecular Weight:** 828.33 g/mol (free acid)**Exact Mass:** 828.09 g/mol (free acid)**Purity:** ≥ 95 % (HPLC)**Form:** solution in water**Color:** orange**Concentration:** 5.0 mM - 5.5 mM**pH:** 7.5 ±0.5**Spectroscopic Properties:** λ<sub>exc</sub> 503 nm, λ<sub>em</sub> 512 nm, ε 80.0 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)**Selected References:**Mancini *et al.* (2007) Exploring the Technology Landscape of 7TMR Drug Signaling Profiling. *Curr. Top. Med. Chem.* **15** (24):2528Vadakkadathmeethal *et al.* (2007) Fluorescence-based adenylyl cyclase assay adaptable to high throughput screening. *Comb. Chem. High Throughput Screen.* **10** (4):289.Kortach *et al.* (2004) Spontaneous nucleotide exchange in low molecular weight GTPases by fluorescently labeled gamma-phosphate-linked GTP analogs. *Proc. Natl. Acad. Sci. USA* **101** (9):2008.Whelan *et al.* (2004) Affinity assays using fluorescence anisotropy with capillary electrophoresis separation. *Anal. Chem.* **76** (24):7380.Gill *et al.* (2003) Low-affinity interactions of BODIPY-FL-GTPγS and BODIPY-FL-GppNHp with G(i)- and G(s)-proteins. *Naunyn Schmiedebergs Arch. Pharmacol.* **368** (3):210.McEwen *et al.* (2001) Fluorescent BODIPY-GTP analogs: real-time measurement of nucleotide binding to G proteins. *Anal. Biochem.* **291** (1):109.