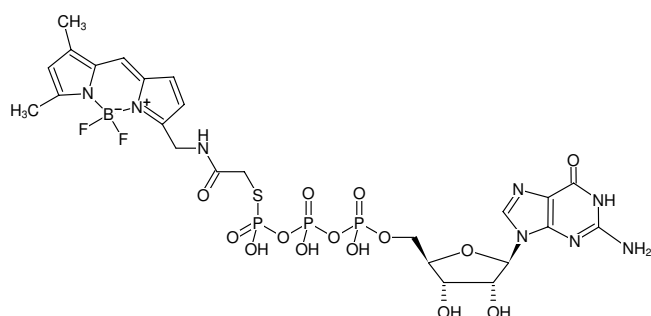


**GTPyS-BDP-FL**

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 gel packs**Storage Conditions:** store at -20 °C**Shelf Life:** 12 months after date of delivery**Molecular Formula:** C₂₄H₃₀BF₂N₈O₁₄P₃S (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_{exc} 503 nm, λ_{em} 512 nm, ε 80.0 L mmol⁻¹ cm⁻¹ (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.