

**ATPyS-BDP-FL**

structural analog to BODIPY® FL

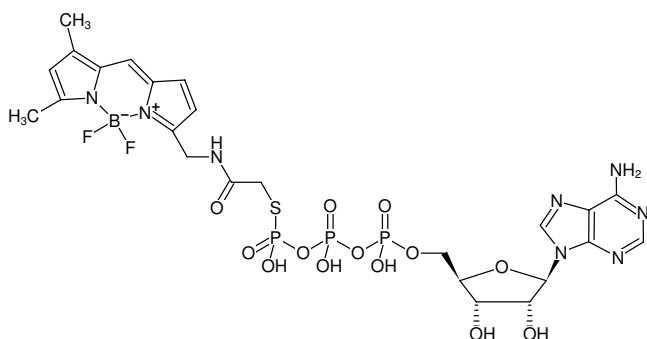
Adenosine-5'-(γ-thio)-triphosphate, labeled with BDP-FL, Sodium salt

Cat. No.	Amount
NU-978	50 µl (5 mM)

Selected References:

Huang *et al.* (2014) Dimeric structure of pseudokinase RNase L bound to 2-5A reveals a basis for interferon-induced antiviral activity. *Mol. Cell.* **53** (2):221.

Kaelyn *et al.* (2012) Activity-Based Probe for Histidine Kinase Signaling *J. Am. Chem. Soc.* **134** (22):9150.



Structural formula of ATPyS-BDP-FL

For general laboratory use.**Shipping:** shipped on gel packs**Storage Conditions:** store at -20 °C**Shelf Life:** 12 months after date of delivery**Molecular Formula:** C₂₄H₃₀BF₂N₈O₁₃P₃S (free acid)**Molecular Weight:** 812.34 g/mol (free acid)**Exact Mass:** 812.09 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)