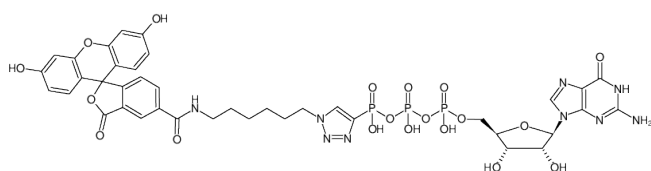




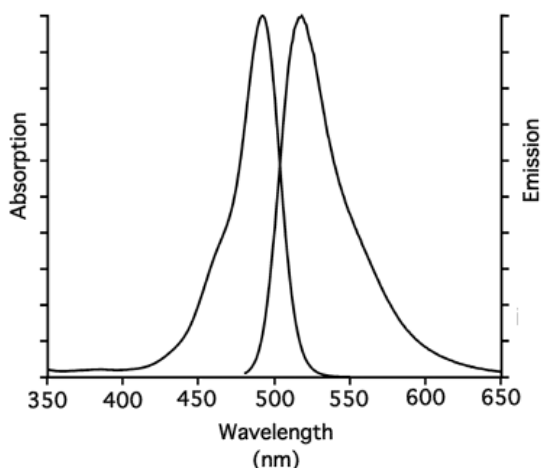
## $\gamma$ -[(6-Aminoethyl)triazolo]-GTP-5-FAM

$\gamma$ -[(6-Aminoethyl)triazolo]- guanosine-5'-triphosphate , labeled with 5-FAM, Triethylammonium salt  
 Guanosine-5'-O-diphosphate[ $\gamma$ -((4-phosphono-1H-1,2,3-triazol-1-yl)hexyl-6-amine-5-FAM)],  
 GTP $\gamma$ -Triazolo-6-AH-5-FAM

| Cat. No.    | Amount            |
|-------------|-------------------|
| NU-1220-5FM | 10 $\mu$ l (1 mM) |



Structural formula of  $\gamma$ -[(6-Aminoethyl)triazolo]-GTP-5-FAM



excitation and emission spectrum of 5-FAM

**For general laboratory use.**

**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>39</sub>H<sub>40</sub>N<sub>9</sub>O<sub>19</sub>P<sub>3</sub> (free acid)

**Molecular Weight:** 1031.71 g/mol (free acid)

**Exact Mass:** 1031.17 g/mol (free acid)

**Purity:**  $\geq$  95 % (HPLC)

**Form:** solution in water

**Color:** yellow to orange

**Concentration:** 1.0 mM - 1.1 mM

**pH:** 7.5  $\pm$  0.5

**Spectroscopic Properties:**  $\lambda_{exc}$  492 nm,  $\lambda_{em}$  517 nm,  
 $\epsilon$  83.0 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)

**Selected References:**

Pepanian *et al.* (2022) A fluorescence anisotropy assay with guanine nucleotides provides access to functional analysis of Gai1 proteins. *Anal Chem.* **94(41)**:14410.