

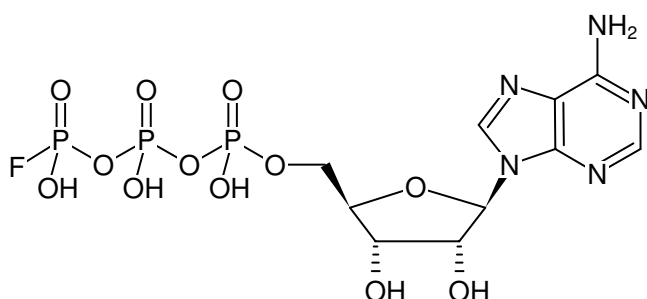
**ATP-y-F**

(ApppF)

Adenosine-5'-(γ-fluoro)-triphosphate, Sodium salt

Adenosine-5'-(3-fluoro)-triphosphate, Sodium salt

Cat. No.	Amount
NU-942-5	5 mg
NU-942-25	25 mg



Structural formula of ATP-y-F

**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>10</sub>H<sub>15</sub>FN<sub>5</sub>O<sub>12</sub>P<sub>3</sub> (free acid)**Molecular Weight:** 509.17 g/mol (free acid)**Exact Mass:** 508.99 g/mol (free acid)**CAS#:** 37515-63-2 (free acid), 1698010-87-5 (sodium salt)**Purity:** ≥ 95 % (HPLC)**Form:** solid**Color:** white to off-white**Spectroscopic Properties:** λ<sub>max</sub> 259 nm, ε 15.4 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)**Applications:**Substrate for snake venom phosphodiesterase <sup>[1,2]</sup>Inhibitor of Lysyl-, Valyl- and Arginyl-tRNA synthases <sup>[3]</sup>Substrate for Fhit proteins <sup>[1,4]</sup>**Selected References:**

[1] Baranowski *et al.* (2015) Synthesis of fluorophosphate nucleotide analogues and their characterization as tools for <sup>19</sup>F NMR studies. *J. Org. Chem.* **80** (8):3982.

[2] Haley *et al.* (1972) g-Fluoroadenosine triphosphate. Synthesis, properties, and interaction with myosin and heavy meromyosin. *Biochemistry* **11** (15):2863.

[3] Freist *et al.* (1980) Chemically modified ATP derivatives for the study of aminoacyl-tRNA synthetases from Bakers' yeast: ATP analogs with fixed conformations of modified triphosphate chains in the aminoacylation reaction *Bioorganic Chemistry* **9** (4):491.

[4] Guranowski *et al.* (2008) Fhit proteins can also recognize substrates other than dinucleoside polyphosphates. *FEBS Lett.* **582** (20):3152.