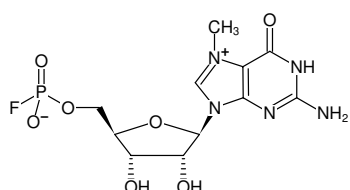


**m<sup>7</sup>GMP-α-F**(m<sup>7</sup>GpF)

7-Methyl-guanosine-5'-(α-fluoro)-monophosphate

7-Methyl-guanosine-5'-(1-fluoro)-monophosphate

Cat. No.	Amount
NU-944-5	5 mg
NU-944-25	25 mg

**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>11</sub>H<sub>15</sub>FN<sub>5</sub>O<sub>7</sub>P**Molecular Weight:** 379.24 g/mol**Exact Mass:** 379.07 g/mol**CAS#:** 1698011-01-6 (inner salt)**Purity:** ≥ 95 % (HPLC)**Form:** solid**Color:** white to off-white**Spectroscopic Properties:** λ<sub>max</sub> 258/280 nm, ε 9.8/8.0 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)**Applications:**

Substrate for human mRNA decapping scavenger enzyme [1,2]

**Selected References:**

[1] Baranowski *et al.* (2016) A fluorescent HTS assay for phosphohydrolases based on nucleoside 5'-fluorophosphates: its application in screening for inhibitors of mRNA decapping scavenger and PDE-I. *Org. Biomol. Chem.* **14** (20):4595.

[2] 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.