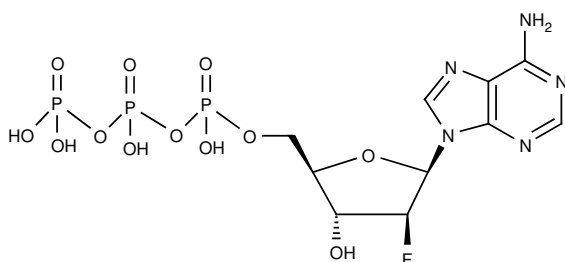


## 2'-Fluoro-2'-deoxy-*ara*-adenine-5'-triphosphate tetra lithium salt

Cat. No.	Amount
NU-10507-A-10	10 $\mu$ mol



**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)

**Storage conditions:** store at -20 °C

**Shelf life:** 12 months

**Purity:** >95 %

**Appearance:** white to off-white powder

**For research use only!**