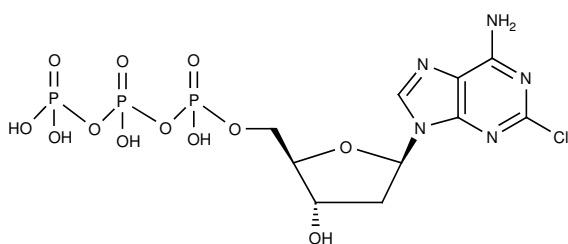


2Cl-dATP, tetra-lithium salt

2'-Deoxy-2-chloroadenosine 5'-triphosphate, tetra-lithium salt

Cat. No.	Amount
NU-10704-10	10 μ mol



Molecular Formula: C₁₀H₁₅ClN₅O₁₂P₃ (free acid)

Molecular Weight: 525,63 g/mol (free acid)

Storage conditions: store at -20 °C

Shelf life: 12 months

Purity: >95 %

Appearance: white to off powder

For research use only!