

Product Information

Description	Fludarabine-Cl has inhibition effect on RNA adenosine deaminase 1 (ADAR1), and can be used for preventing and/or researching cancer or tumor-related diseases.
Molecular Formula:	C ₁₀ H ₁₁ ClFN ₅ O ₃
Supplied As:	White powder
Solubility:	H ₂ O : 5 mg/ml (16.46 mM; ultrasonic and warming and heat to 60°C)
Stability:	Store solid as supplied at -20°C for up to 3 years. Store desiccated as supplied at +4°C for up to 2 years. Store solutions at -80°C for up to 6 months. Store solutions at -20°C for up to 1 month. Avoid freeze/thaw cycles.
Storage:	-20°C
MW:	303.68 Da
Cas No.	2734853-80-4
Purity:	≥99%

Chemical Structure

