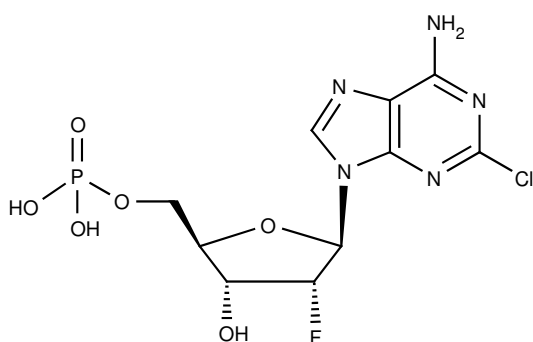


Clofarabine-5'-monophosphate Triethylammonium salt

Cat. No.	Amount
NU-OXEL007-2	3 mg



Molecular Formula: C₁₀H₁₂ClFN₅O₆P

Molecular Weight: 383.66 g/mol (free acid)

Storage conditions: store at -20°C

Shelf life: 12 months

Purity: >95% (HPLC)

Appearance: white solid

For research use only!

Selected References:

Nagai *et al.* (2011) Deoxycytidine kinase modulates the impact of the ABC transporter ABCG2 on clofarabine cytotoxicity. *Cancer Res* **71(5)**:1781.

Xie *et al.* (1995) Metabolism and actions of 2-chloro-9-(2-deoxy-2-fluoro-β-D-arabinofuranosyl)adenine in human lymphoblastoid cells. *Cancer Research* **55(13)**:2847.