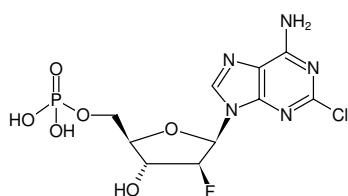


**Clofarabine-5'-monophosphate**

Triethylammonium salt

| Cat. No. | Amount |
|----------|--------|
| NU-873   | 3 mg   |



Structural formula of Clofarabine-5'-monophosphate

**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>10</sub>H<sub>12</sub>ClFN<sub>5</sub>O<sub>6</sub>P (free acid)**Molecular Weight:** 383.66 g/mol (free acid)**Exact Mass:** 383.02 g/mol (free acid)**CAS#:** 134646-42-7**Purity:** ≥ 95 % (HPLC)**Form:** solid**Color:** white to off-white**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.