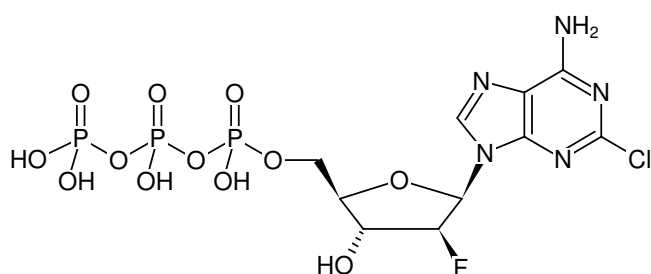




## Clofarabine-5'-triphosphate

Triethylammonium salt

Cat. No.	Amount
NU-874	5 $\mu$ mol



Structural formula of Clofarabine-5'-triphosphate

**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>14</sub>ClFN<sub>5</sub>O<sub>12</sub>P<sub>3</sub> (free acid)

**Molecular Weight:** 543.62 g/mol (free acid)

**Exact Mass:** 542.95 g/mol (free acid)

**CAS#:** 167620-89-5

**Purity:**  $\geq$  95 % (HPLC)

**Form:** oil

**Color:** colorless to slightly yellow

**Spectroscopic Properties:**  $\lambda_{\max}$  262 nm,  $\epsilon$  14.3 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)

### Selected References:

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

Yamauchi *et al.* (2001) DNA repair initiated in chronic lymphocytic leukemia lymphocytes by 4-hydroperoxycyclophosphamide is inhibited by fludarabine and clofarabine. *Clinical Cancer Research* **7** (11):3580.

Parker *et al.* (1991) Effects of 2-chloro-9-(2-deoxy-2-fluoro- $\beta$ -D-arabinofuranosyl)adenine on K562 cellular metabolism and the inhibition of human ribonucleotide reductase and DNA polymerases by its 5'-triphosphate. *Cancer Research* **51** (9):2386.