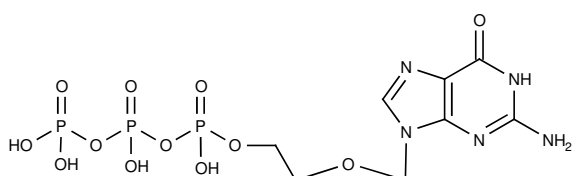


## Aciclovir-5'-triphosphate Triethylammonium salt

Cat. No.	Amount
NU-OXEL020-2	2 mg



**CAS number:** 66341-18-2

**Molecular Formula:** C<sub>8</sub>H<sub>14</sub>N<sub>5</sub>O<sub>12</sub>P<sub>3</sub>

**Molecular Weight:** 465.14 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:

Keller *et al.* (2008) Interaction of Herpes Primase with the Sugar of a NTP. *Biochemistry* **47**(34):8977.

Hanes *et al.* (2007) Enzymatic Therapeutic Index of Acyclovir: Viral versus human polymerase  $\gamma$  specificity. *Journal of Biological Chemistry* **282**(34):25159.

Huang *et al.* (1999) The enzymological basis for resistance of herpesvirus DNA polymerase mutants to acyclovir: relationship to the structure of  $\alpha$ -like DNA polymerases. *Proceedings of the National Academy of Sciences of the United States of America* **96**(2):447.