

**dITP - Solid**

2'-Deoxyinosine-5'-triphosphate, Sodium salt

Cat. No.	Amount
NU-1040-20	20 µmol
NU-1040-100	5 x 20 µmol

**For research use only!****Shipping:** shipped on blue ice**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>15</sub>N<sub>4</sub>O<sub>13</sub>P<sub>3</sub> (free acid)**Molecular Weight:** 492.16 g/mol (free acid)**CAS#:** 16595-02-1**Purity:** ≥ 95 % (HPLC)**Form:** lyophilised**Spectroscopic Properties:** λ<sub>max</sub> 249 nm, ε 12.2 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.0)**Selected References:**Erlich *et al.* (1988) Primer-directed enzymatic amplification of DNA with a thermostable DNA polymerase. *Science* **29** (239):487.Holland *et al.* (1991) Detection of specific polymerase chain reaction product by utilizing the 5'→3' exonuclease activity of *Thermus aquaticus* DNA polymerase. *Proc. Natl. Acad. Sci. USA* **88** (16):7276.Sanger *et al.* (1977) DNA sequencing with chain-terminating inhibitors. *Proc. Natl. Acad. Sci. USA* **74**:5463.