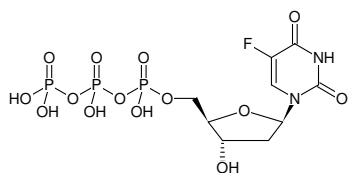


**5-Fluoro-dUTP**

(5F-dUTP)

5-Fluoro-2'-deoxyuridine-5'-triphosphate, Sodium salt

Cat. No.	Amount
NU-154S	20 µl (100 mM)
NU-154L	5 x 20 µl (100 mM)

**Selected References:**Longley *et al.* (2003) 5-Fluorouracil: mechanisms of action and clinical strategies. *Nature Reviews Cancer*. **3**:300.

Structural formula of 5-Fluoro-dUTP

**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>9</sub>H<sub>14</sub>N<sub>2</sub>O<sub>14</sub>P<sub>3</sub>F (free acid)**Molecular Weight:** 486.13 g/mol (free acid)**Exact Mass:** 485.96 g/mol (free acid)**Purity:** ≥ 95 % (HPLC)**Form:** clear aqueous solution**Concentration:** 100 mM - 110 mM**pH:** 7.5 ± 0.5**Spectroscopic Properties:** λ<sub>abs</sub> 267 nm, ε 10.0 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)