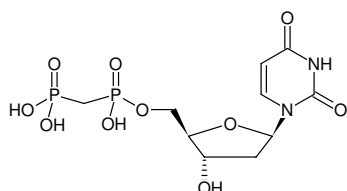


**dUpCp**

(dUMPCP)

2'-Deoxyuridine-5'-[( $\alpha,\beta$ )-methylene]diphosphate, Sodium salt

Cat. No.	Amount
NU-428S	100 $\mu$ l (10 mM)
NU-428L	5 x 100 $\mu$ l (10 mM)



Structural formula of dUpCp

**For research use only!****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>16</sub>N<sub>2</sub>O<sub>10</sub>P<sub>2</sub> (free acid)**Molecular Weight:** 386.19 g/mol (free acid)**Exact Mass:** 386.03 g/mol (free acid)**Purity:**  $\geq$  95 % (HPLC)**Form:** solution in water**Color:** colorless to slightly yellow**Concentration:** 10 mM - 11 mM**pH:** 7.5  $\pm$  0.5**Spectroscopic Properties:**  $\lambda_{\max}$  262 nm,  $\epsilon$  10.2 L mmol<sup>-1</sup> cm<sup>-1</sup>  
(Tris-HCl pH 7.5)