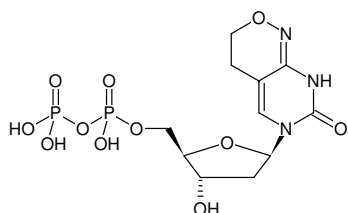


**dPDP**

6H,8H-3,4-Dihydro-pyrimido(4,5-c)(1,2)oxazin-7-one-8-β-D-2'-deoxy-ribofuranoside-5'-diphosphate,
Sodium salt

Cat. No.	Amount
NU-1159S	20 µl (10 mM)
NU-1159L	5 x 20 µl (10 mM)



Structural formula of dPDP

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: C₁₁H₁₇N₃O₁₁P₂ (free acid)

Molecular Weight: 429.21 g/mol (free acid)

Exact Mass: 429.03 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

Form: solution in water

Color: colorless to slightly yellow

Concentration: 10 mM - 11 mM

pH: 7.5 ±0.5