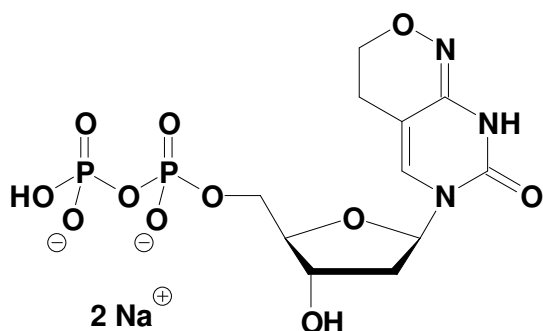


dPDP

6H,8H-3,4-Dihydro-Pyrimido[4,5-c][1,2]oxazin-7-one-8- β -D-2'-Deoxy-Ribofuranosid-5'-diphosphate, Sodium salt

Cat. No.	Amount
NU-1159S	20 Units
NU-1159L	100 Units



Cat. No.: NU-1159

Molecular Formula: C₁₁H₁₅N₃O₁₁P₂ (Anion)

Molecular Weight: 427.20 (Anion)

Purity: > 95%, clear aqueous solution, pH 7.5

Storage conditions:

Short term exposure (up to 1 week cumulative) to ambient temperature possible. Long term storage at < -20°C. If stored as recommended, Jena Bioscience guarantees optimal performance of this product for 12 months after date of delivery.

* 1 unit = 1 μ l of a 10 mM solution

For research use only!