

# Sp-dPTP- $\alpha$ -S

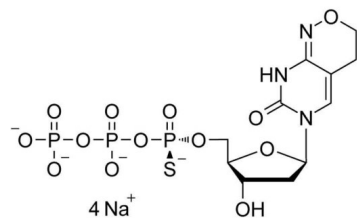
Triphosphate analogue with artificial base for directed mutagenesis experiments. Increased metabolic stability compared to dPTP (Cat. No. D 114). Accepted by DNA polymerase for incorporation of phosphorothioate into DNA.

*BLG-D106-05 (5 x 1  $\mu$ mol pack size) is not sold in the U.S. or Canada.  
Please [contact us](#) for available options.*

## Ordering Information [Order Online »](#)

BLG-D106-01	1 $\mu$ mol
BLG-D106-05	5x1 $\mu$ mol

## Manuals, SDS & CofA [View Online »](#)



## Handling & Storage

Long Term Storage -80°C

Shipping Dry Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** 6-(β-D-2-Deoxyribofuranosyl)-3, 4-dihydro-8H-pyrimido-[4, 5-c]-[1, 2]-oxazin-7-one- 5'-O-(1-thiotriphosphate) . Sp- isomer, 2'-Deoxy-P-nucleoside-5'-O-(1-thiotriphosphate) . Sp-isomer

**CAS** 627502-68-5

**Couple Type** Modified nucleotides

**Formula**  $C_{11}H_{18}N_3O_{13}P_3S$  (free acid)

**Formulation** 10 mM aqueous solution of pH 7.6

**MW** 525.3

**Purity** > 95% HPLC

**Quantity** 1 μmol is equivalent to ~0.5 mg



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-  
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-  
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-  
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-  
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-  
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)