

pdCp is accepted by T4 RNA ligase to build oligodeoxyribonucleotides.

*BLG-D094-25 (5 x 5  $\mu$ mol pack size) is not sold in the U.S. or Canada.*

Please [contact us](#) for available options.

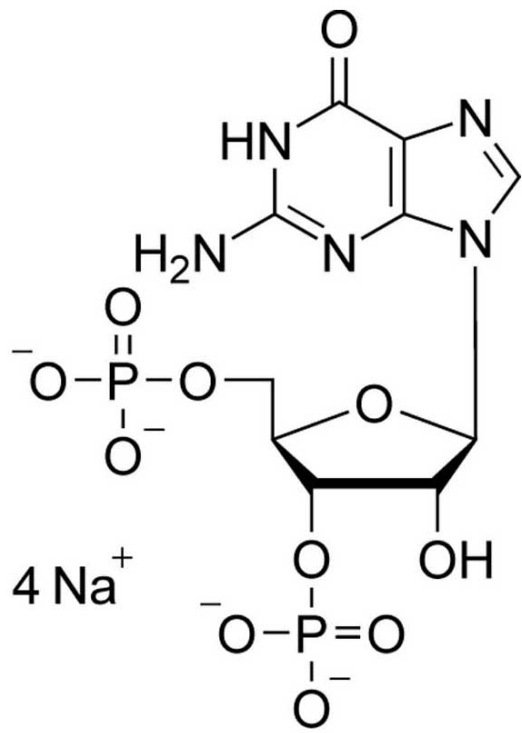
## Ordering Information

[Order Online »](#)

|             |               |
|-------------|---------------|
| BLG-D094-05 | 5 $\mu$ mol   |
| BLG-D094-25 | 5x5 $\mu$ mol |

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

Long Term Storage -20°C

Shipping Ambient Temperature

**Regulatory Status** RUO - Research Use Only

## Product Details

Alternative Name 2'-Deoxyguanosine-3',5'-O-bisphosphate . sodium salt

CAS 16174-59-7

Couple Type Modified nucleotides

Formula  $C_{10}H_{15}N_5O_{10}P_2$  (free acid)

MW 427.2 (free acid)

Purity > 95% HPLC

Quantity 5µmol≈2.1mg

Solubility Soluble in water (≥36mM).

**Technical Info / Product Notes** For the Original Manufacturer's data sheet please [click here](#).



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)