

# 2'-/3'-AHC-GDP

Functionalized GDP analogue for immobilization or conjugation with markers and dyes

2'-/3'-AHC-GDP is an analogue of guanosine-5'-O-diphosphate (GDP) in which a hexyl spacer with a terminal amino group has been attached either to the ribose 2'-hydroxy group or to the 3'-hydroxy group by a carbamate bond. It is suitable as a ligand in affinity chromatography or for coupling of various labelling structures including fluorophores. The spacer is connected to the ribose 2'-hydroxy group of the guanosine part.

*BLG-A289-25 (5 x 5 µmol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.*

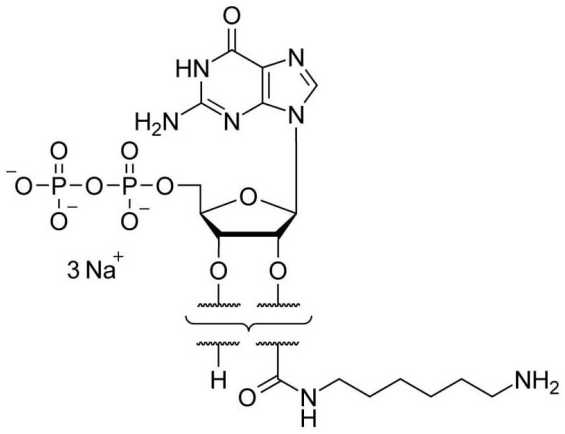
## Ordering Information

[Order Online »](#)

BLG-A289-05	5µmol
BLG-A289-25	5x5µmol

## 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** 2'-/3'-O-(6-Aminohexylcarbamoyl)guanosine-5'-O-diphosphate . sodium salt

**CAS** 1810735-99-9 / 1810737-23-5

**Formula**  $C_{17}H_{29}N_7O_{12}P_2$  (free acid)

**Formulation** 10 mM aqueous solution

**MW** 585.4 (free acid)

**Purity** > 95% HPLC

**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)