

# 6-T-GDP

6-T-GDP is a metabolite of azathioprine, an immunosuppressive drug that is used in the treatment of several diseases, including childhood acute lymphoblastic leukaemia (ALL) and inflammatory bowel disease.

*BLG-T015-25 (5 x 5 µmol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.*

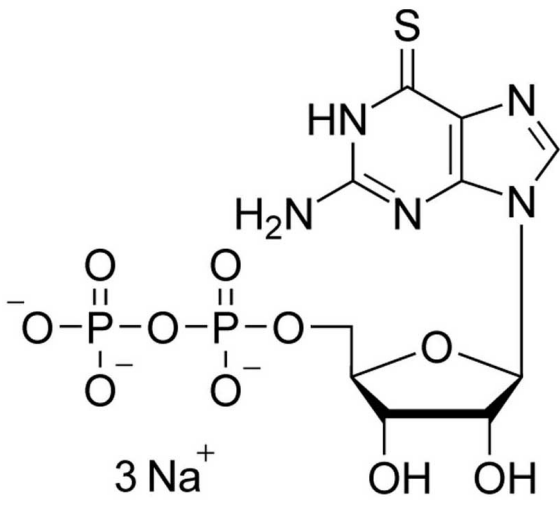
## Ordering Information

[Order Online »](#)

BLG-T015-05	5µmol
BLG-T015-25	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** 6-Thioguanosine-5'-O-diphosphate . trisodium salt

**Appearance** 10 mM aqueous solution

**CAS** 16541-19-8

**Couple Type** Modified nucleotides

**Formula**  $C_{10}H_{15}N_5O_{10}P_2S$  (free acid)

**MW** 459.3 (free acid)

**Purity** > 95% HPLC

**Quantity** 5µmol≈2.3mg

**Solubility** Soluble in water.

**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)