

# 5'-dIMPS

Potential substrate, competitive inhibitor or regulator of enzymes that interacts with 2'-deoxyinosine-5'-monophosphate.

*BLG-D060-25 (5 x 5 µmol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.*

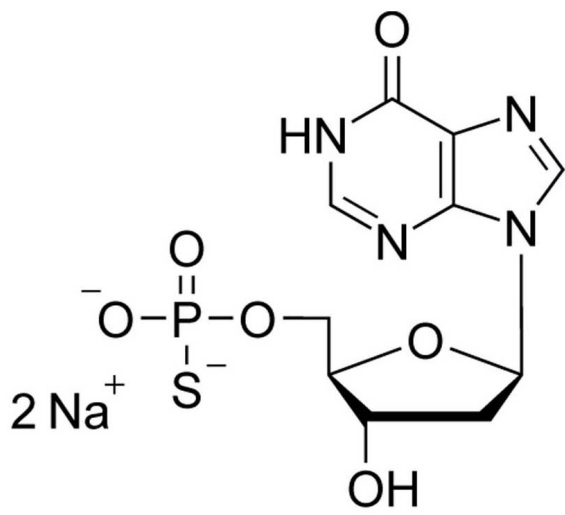
## Ordering Information

[Order Online »](#)

BLG-D060-05	5µmol
BLG-D060-25	5x5µmol

## 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'-Deoxyinosine-5'-O-monophosphorothioate . disodium salt

CAS 771477-45-3

Couple Type Modified nucleotides

Formula  $C_{10}H_{13}N_4O_6PS$  (free acid)

MW 348.3 (free acid)

Purity > 97% HPLC

Quantity 5µmol≈1.7mg



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)