

# CI-ANT-dATP

## Fluorescent dATP analogue

Analogue of 2'-deoxyadenosine-5'-triphosphate with only moderate intrinsic fluorescence ( $\lambda_{exc}$  347 nm,  $\lambda_{em}$  ~431 nm). The fluorophore is supposed to have a certain sensitivity for its environment and can change its spectral properties upon binding. Useful for research into dATP-dependent receptor proteins. Potential inhibitor of various adenylyl cyclase isoforms and soluble guanylyl cyclase (sGC).

*BLG-C142-05 (5 x 1  $\mu$ mol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.*

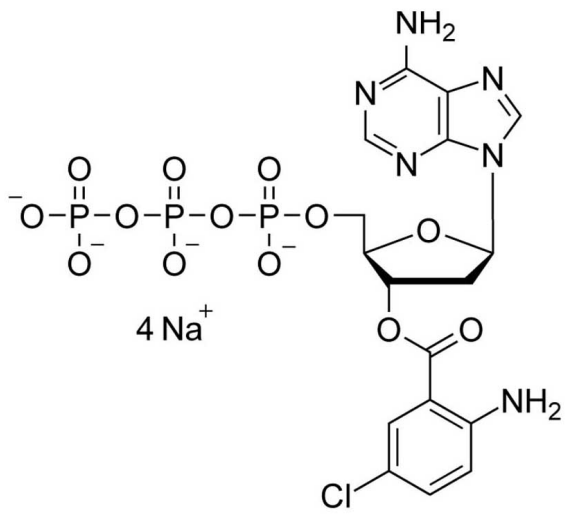
### Ordering Information

[Order Online »](#)

BLG-C142-01	1 $\mu$ mol
BLG-C142-05	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** 3'-O-(5-Chloroanthraniloyl)-2'-deoxyadenosine-5'-O-triphosphate . sodium salt

**Couple Type** Modified nucleotides

**Formula**  $C_{17}H_{20}ClN_6O_{13}P_3$  (free acid)

**Formulation** 10 mM aqueous solution

**MW** 644.8 (free acid)

**Purity** > 95% HPLC

**Quantity** 1  $\mu$ mol is equivalent to ~0.6 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)