

# CI-ANT-ATP

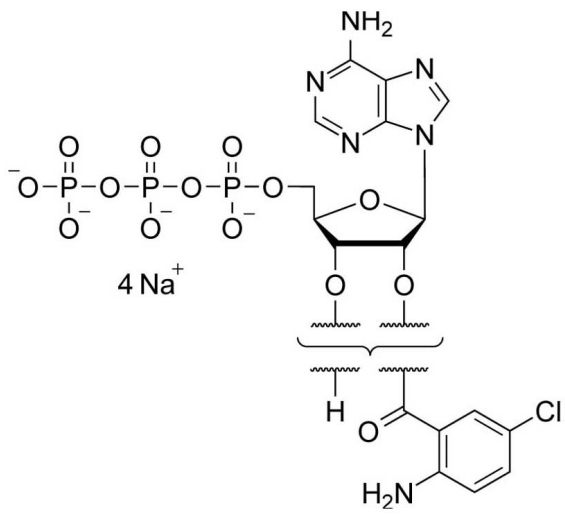
Analogue of adenosine-5'-triphosphate with only moderate intrinsic fluorescence ( $\lambda_{exc}$  350 nm;  $\lambda_{em}$  ~450 nm), increasing considerably within hydrophobic environment. Useful for research into ATP-dependent receptor proteins.

*BLG-C114-25 (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-C114-05	5 $\mu$ mol
BLG-C114-25	5x5 $\mu$ mol

## Manuals, SDS & CofA [View Online »](#)



## Handling & Storage

**Handling** Protect from light.

**Long Term Storage** -80°C

**Shipping** Dry Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** 2'-/3'-O-(5-Chloroanthraniloyl)adenosine-5'-O-triphosphate . tetrasodium salt

**Appearance** 10 mM aqueous solution

**CAS** 1350521-45-7

**Couple Type** Modified nucleotides

**Formula**  $C_{17}H_{20}ClN_6O_{14}P_3$  (free acid)

**MW** 660.8

**Purity** > 95% HPLC



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)