

# 8-N<sub>3</sub>-ε-ATP

8-N<sub>3</sub>-ε-ATP combines the photosensitive activity of 8-N<sub>3</sub>-ATP (Cat. No. A 043) and the fluorescent properties of ε-ATP (Cat. No. E 004). It is useful for photoaffinity labelling of ATP binding proteins and subsequent detection of the fluorescent label ( $\lambda_{exc}$  285 nm,  $\lambda_{em}$  412 nm). Alternatively, detection may be achievable by using an ethenoadenosine-specific antibody (1G4, compare: Krebs et al., Anal. Biochem., 314, 108 - 115 (2003)).

*BLG-A211-25 (5 x 5  $\mu$ mol pack size) is not sold in the U.S. or Canada.  
Please [contact us](#) for available options.*

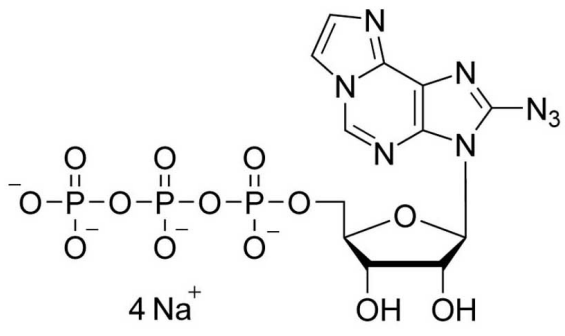
## Ordering Information

[Order Online »](#)

BLG-A211-05	5 $\mu$ mol
BLG-A211-25	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** 8-Azido-1, N6-ethenoadenosine-5'-O-triphosphate . tetrasodium salt

**Appearance** 10 mM aqueous solution

**CAS** 66895-05-4

**Couple Type** Modified nucleotides

**Formula**  $C_{12}H_{15}N_8O_{13}P_3$  (free acid)

**MW** 572.2

**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)