

# Etheno-ADP ( $\epsilon$ -ADP), L pack

## Ordering Information

[Order Online »](#)

JBS-NU-1140L	5x500 $\mu$ l
--------------	---------------

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** 1,N6-Etheno-ADP, 1,N6-Etheno-adenosine-5'-diphosphate, Sodium salt

**Appearance** colorless to slightly yellow solution

**CAS** 38806-39-2

**Couple Type** Modified nucleotides

**Formula**  $C_{12}H_{15}N_5O_{10}P_2$  (free acid)

**Formulation** solution in water

**MW** 451.22 (free acid)

**Purity** ≥ 95 % (HPLC)

**Quantity** 1 unit = 1 µl of a 10 mM solution

**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)