

# EDA-ADP – ATTO-MB2

## Ordering Information

[Order Online »](#)

|                |      |
|----------------|------|
| JBS-NU-802-MB2 | 40µl |
|----------------|------|

## 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                    2’/3’-O-(2-Aminoethyl-carbamoyl)-Adenosine-5’-diphosphate, labeled with ATTO-MB2, Triethylammonium salt

Couple Type                        Modified nucleotides

Formula                              C<sub>32</sub>H<sub>40</sub>N<sub>10</sub>O<sub>12</sub>P<sub>2</sub>S (free acid)

Formulation                        solution in water

MW                                    850.73 (free acid)

Purity                                ≥ 95 % (HPLC)

Technical Info / Product Notes                    For the Original Manufacturer’s data sheet please [click here](#).

Image not found or type unknown  
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)