

# Adenosine 5'- monophosphate . disodium salt

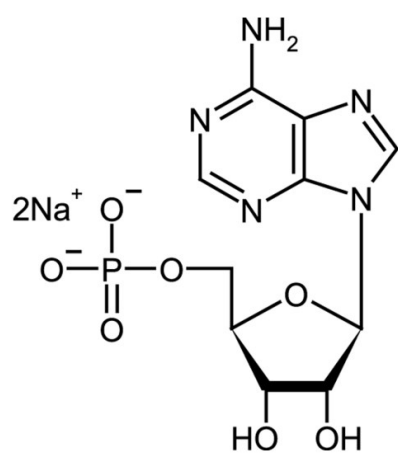
## Ordering Information

[Order Online »](#)

|                  |    |
|------------------|----|
| ALX-480-014-G005 | 5g |
|------------------|----|

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

|                   |                                                                                                                                                                                                              |
|-------------------|--------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------|
| Use/Stability     | As indicated on product label or CoA when stored as recommended. Solutions should be stored in the refrigerator and should be lyophilized and frozen for longer storage periods. The product is not sterile. |
| Handling          | Protect from moisture.                                                                                                                                                                                       |
| Long Term Storage | -20°C                                                                                                                                                                                                        |
| Shipping          | Ambient Temperature                                                                                                                                                                                          |

## Regulatory Status

RUO - Research Use Only

## Product Details

|                  |                                      |
|------------------|--------------------------------------|
| Alternative Name | AMP . 2Na                            |
| Appearance       | White to off-white powder.           |
| CAS              | 4578-31-8                            |
| Formula          | $C_{10}H_{12}N_5O_7P \cdot 2Na$      |
| MI               | 14: 158                              |
| MW               | 345.2 . 46.0                         |
| Purity           | ≥98% (HPLC)                          |
| RTECS            | AU7481600                            |
| Solubility       | Soluble in water or aqueous buffers. |

Last modified: May 29, 2024



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-  
fr@enzolifesciences.com](mailto:info-fr@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)