



Control reagent for nucleotide-AM ester applications

Membrane-permeable prodrug of inorganic phosphate. The acetoxymethyl groups mask the charged polar phosphate and thus render the molecule highly membrane-permeant. Inside the cell esterases release the highly polar inorganic phosphate (PO_4^{3-}). In addition, the AM ester groups liberate acetic acid and formaldehyde, two metabolites with potential biological functions.

$\text{PO}_4\text{-AM}_3$ is recommended as control reagent in nucleotide-AM ester applications to test for side effects of acetic acid and formaldehyde.

BLG-P030-015 (5 x 0.33 μmol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.

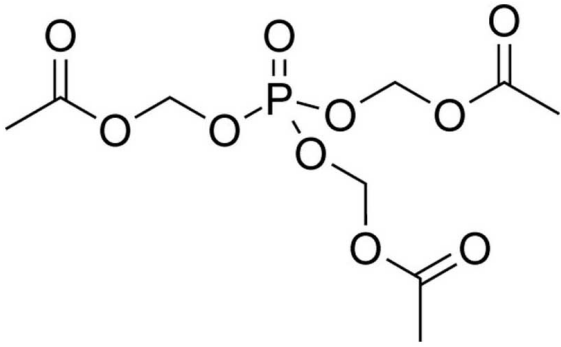
Ordering Information

[Order Online »](#)

BLG-P030-015	5x0.33 μmol
--------------	------------------------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability PO4-AM3 is rather labile in aqueous solution. Therefore, aqueous solutions should be freshly prepared immediately before use. Stock solutions in anhydrous DMSO or DMF should be relatively stable when stored frozen at -20°C to -70°C.

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Phosphate tris(acetoxymethyl)ester

CAS 148334-61-6

Couple Type Substrate

Formula $C_9H_{15}O_{10}P$

MW 314.2

Purity > 95% (MS)

Solubility Due to its rather high lipophilicity, the solubility of PO4-AM3 in water or buffers is limited. We suggest to use a small amount of anhydrous organic solvent such as anhydrous DMSO or DMF for preparation of 1-100mM stock solutions, and to dilute with water or buffer down to the concentrations required. In some cases, especially at high concentrations (~ 1mM), Pluronic® F-127 (Molecular Probes) can be useful to facilitate solubilization in water or physiological media. Please be sure to check for DMSO/DMF tolerance in your system. Since PO4-AM3 is bioactivated by esterases, application to cell cultures should be performed without serum supplements (even heat-inactivated serum still contains active esterases!) in the media for at least 15 minutes. Otherwise, serum esterases may strongly reduce the cell-loading efficacy. Rinse tube walls carefully and preferably use ultrasonic or vortex to achieve total and uniform mixing.

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)