

# Ap4A /AppppA

Induces proliferation of vascular smooth muscle cells (Bobbert et al., Biochem. Pharmacol., 75, 1966 - 1973 (2008)), apoptosis in human cultured cells (Vartanian et al., FEBS Lett., 456, 175 - 180 (1999)) and exhibits vasoconstrictor activity (Lewis et al., Br. J. Pharmacol., 129, 124 - 130 (2000)).

*BLG-D115-05 (5 x 1  $\mu$ mol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.*

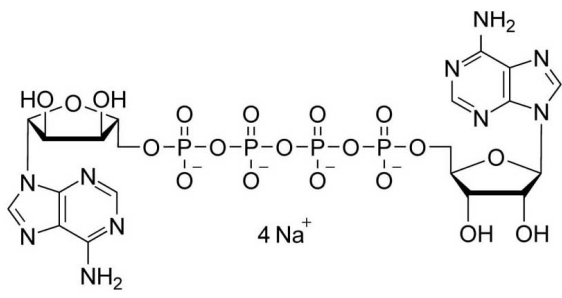
## Ordering Information

[Order Online »](#)

BLG-D115-01	1 $\mu$ mol
BLG-D115-05	5x1 $\mu$ mol

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

Long Term Storage -20°C

Shipping Ambient Temperature

**Regulatory Status** RUO - Research Use Only

## Product Details

Alternative Name P1, P4-Di-(adenosine-5')-tetraphosphate . tetrasodium salt

CAS 5542-28-9

Couple Type Modified nucleotides

Formula  $C_{20}H_{28}N_{10}O_{19}P_4$  (free acid)

MW 836.4

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)