

2'/3'-Biotin-16-Ap₄A

Biotinylated analogue of Ap₄A

BLG-B289-005 (5 x 0.1 µmol pack size) is not sold in the U.S. or Canada.

Please [contact us](#) for available options.

Ordering Information

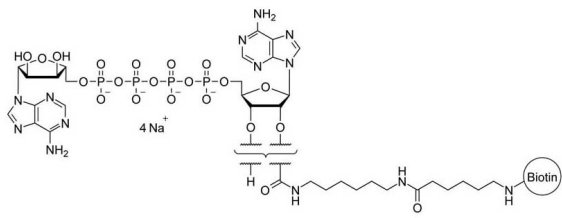
[Order Online »](#)

BLG-B289-001	0.1 µmol
BLG-B289-005	5x0.1 µmol

Manuals, SDS & CofA

[View Online »](#)

Biotinylated analogue of the second messenger P¹,P⁴-Di-(adenosine-5')-tetraphosphate (Ap₄A, Cat. No. D 115). Potential tracer in immunoassays, for affinity chromatography or pull-down assays.



Handling & Storage

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name P1-(5'-[2'-/ 3'-O-(6-(6-[Biotinyl]aminohexanoyl)aminohexylcarbamoyl)adenosyl])-P4-(5'-adenosyl)-tetraphosphate . sodium salt, 2'/3'-Biotin-16-AppppA

Excitation Maximum 259 nm

Formula $C_{43}H_{67}N_{15}O_{23}P_4S$ (free acid)

MW 1318.1 (free acid)

Purity > 95% HPLC

Quantity 0.1 µmol / ~ 132 µg



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com