

ϵ -Ap₄- ϵ -A

Fluorescent analogue of Ap₄A

Fluorescent analogue of the parent compound Ap₄A / AppppA (Cat. No. BLG-D115) with λ_{exc} 307 nm, λ_{em} 410 nm.

BLG-E030-05 (5 x 1 μ mol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.

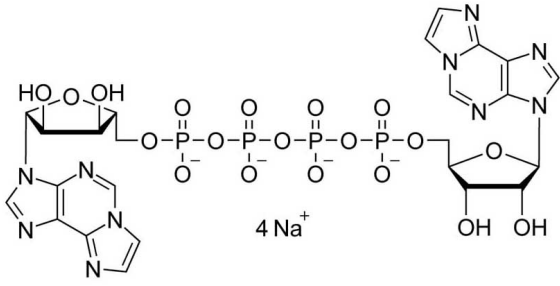
Ordering Information

[Order Online »](#)

| | |
|-------------|---------------|
| BLG-E030-01 | 1 μ mol |
| BLG-E030-05 | 5x1 μ 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 ϵ -Apppp- ϵ -A, P1-(5'-[1,N6-Ethenoadenosyl])-P4-(5'-[1,N6-Ethenoadenosyl])-tetraphosphate . sodium salt

Appearance Solid

CAS 96920-51-3

Formula $C_{24}H_{28}N_{10}O_{19}P_4 \cdot 4 Na$

MW 884.4

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

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