

ϵ -Ap₄U

Fluorescent analogue of Up₄A

Fluorescent analogue of the purinoceptor agonist Up₄A (Prod. No. BLG-U008); λ_{exc} 230 nm, λ_{em} 420 nm.

*BLG-E031-05 (5 x 1 μ mol pack size) is not sold in the U.S. or Canada.
Please [contact us](#) for available options.*

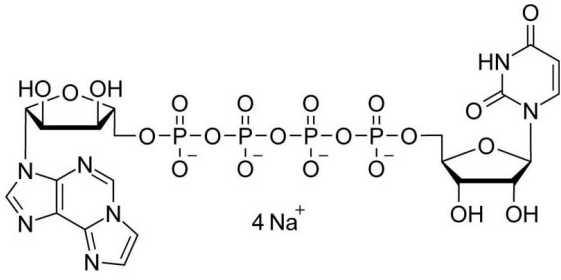
Ordering Information

[Order Online »](#)

BLG-E031-01	1 μ mol
BLG-E031-05	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 -AppppU, P1-(5'-[1, N6-Ethenoadenosyl])-P4-(5'-uridy)-tetraphosphate, ϵ -Up4A

Formula $C_{21}H_{27}N_7O_{21}P_4$ (free acid)

MW 837.4 (free acid)

Purity > 95% HPLC



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