

ϵ -5'-AMP

Fluorescent analogue of 5'-AMP

Fluorescent analogue of adenosine-5'-monophosphate; λ_{exc} 300 nm, λ_{em} 415 nm.

*BLG-E005-50 (5 x 10 μ mol pack size) is not sold in the U.S. or Canada.
Please [contact us](#) for available options.*

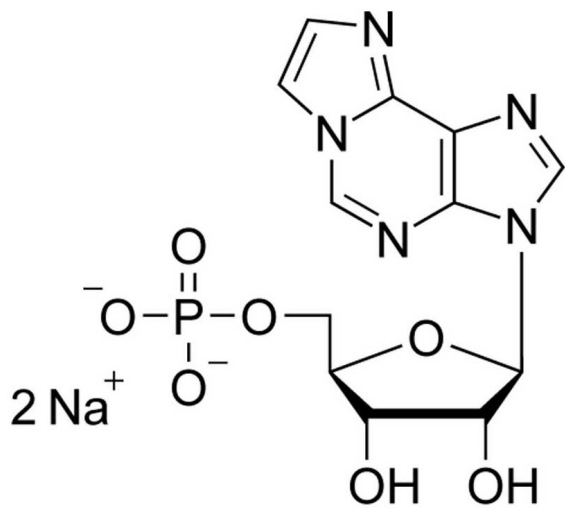
Ordering Information

[Order Online »](#)

BLG-E005-50	5x10 μ mol
-------------	----------------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Handling	Protect from light.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name 1,N6-Ethenoadenosine-5'-O-monophosphate . disodium salt

CAS 103213-41-8

Couple Type Modified nucleotides

Formula $C_{12}H_{14}N_5O_7P$ (free acid)

MW 371.3

Purity > 97% HPLC

Quantity 50µmol≈~21mg

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