

ε-ADPR

Fluorescent ADPR analogue

ε-ADPR is a fluorescent analogue of ADP-ribose ($\lambda_{\text{ex}}=300$ nm and $\lambda_{\text{em}}=410$ nm) and can be used in fluorometric assays.

BLG-E013-50 (5 x 10 µmol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.

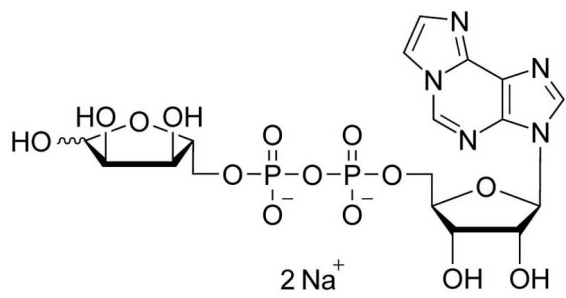
Ordering Information

[Order Online »](#)

BLG-E013-10	10µmol
BLG-E013-50	5x10µmol

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-diphosphoribose . sodium salt
CAS	69699-76-9
Couple Type	Modified nucleotides
Formula	$C_{17}H_{23}N_5O_{14}P_2$ (free acid)
MW	583.3 (free acid)
Purity	> 95% HPLC
Quantity	10µmol≈~5.8mg
Solubility	Soluble in water (≥20mM).

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