

ϵ -5'-dAMP

Fluorescent analogue of 2'-deoxy-5'-AMP

Fluorescent analogue of 2'-deoxyadenosine-5'-monophosphate; λ_{exc} 300 nm, λ_{em} 415 nm.

*BLG-D077-25 (5 x 5 μ mol pack size) is not sold in the U.S. or Canada.
Please [contact us](#) for available options.*

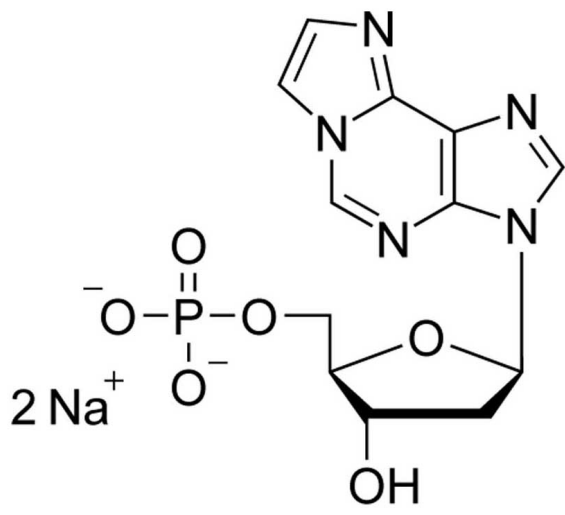
Ordering Information

[Order Online »](#)

BLG-D077-25	5x5 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	2'-Deoxy-1, N6-ethenoadenosine-5'-O-monophosphate . sodium salt
CAS	60508-81-8
Couple Type	Modified nucleotides
Formula	$C_{12}H_{14}N_5O_6P$ (free acid)
MW	355.3
Purity	> 97% HPLC
Quantity	5µmol≈1.8mg
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-fr@enzolifesciences.com

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

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

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