

8-N₃-ε-ADP

8-N₃-ε-ADP combines the photosensitive activity of 8-N₃-ADP (Cat. No. A 044) and the fluorescent properties of ε-ADP (Cat. No. E 007). It is useful for photoaffinity labelling of ADP binding proteins and subsequent detection of the fluorescent label (λ_{exc} 285 nm, λ_{em} 410 nm). Alternatively, detection may be achievable by using an ethenoadenosine-specific antibody (1G4, compare: Krebs et al., Anal. Biochem., 314, 108 - 115 (2003))

*BLG-A210-25 (5 x 5 μ mol pack size) is not sold in the U.S. or Canada.
Please [contact us](#) for available options.*

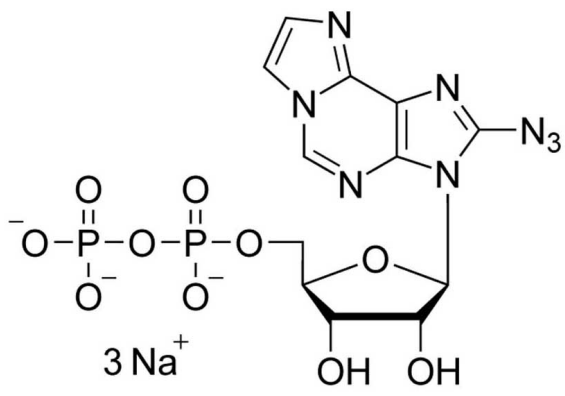
Ordering Information

[Order Online »](#)

BLG-A210-05	5 μ mol
BLG-A210-25	5x5 μ mol

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Handling Protect from light.

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name 8-Azido-1, N6-ethenoadenosine-5'-O-diphosphate . trisodium salt

Appearance 10 mM aqueous solution

Couple Type Modified nucleotides

Formula $C_{12}H_{15}N_8O_{10}P_2$ (free acid)

MW 493.3

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