

6-Fluo-10-NAD⁺

Fluorescent analogue of NAD⁺

6-Fluo-10-NAD⁺ is an analogue of β -NAD⁺ in which 5(6)-carboxyfluorescein has been attached to the amino group in position 6 of the adenine nucleobase via a 10-atom spacer. 6-Fluo-10-NAD⁺ is a fluorescent analogue of β -NAD⁺ (λ_{exc} 494 nm, λ_{em} 517 nm) which can provide a non-radioactive alternative to radio-labelled β -NAD⁺ for research on enzymes requiring β -NAD⁺ as a substrate or cofactor.

BLG-N023-005 (5 x 0.1 μ mol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.

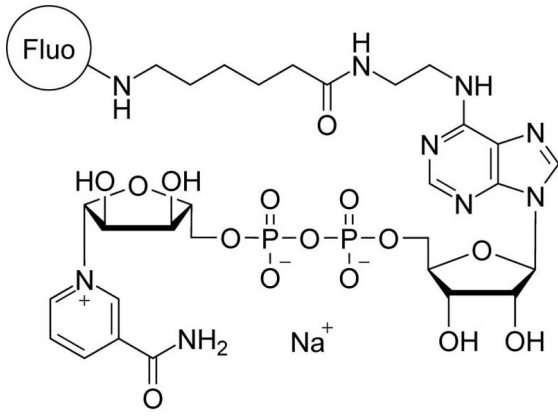
Ordering Information

[Order Online »](#)

BLG-N023-001	0.1 μ mol
BLG-N023-005	5x0.1 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 β -Nicotinamide-N6-(2-(6-[fluoresceinyl]aminohexanoyl)aminoethyl)adenine dinucleotide . sodium salt

Couple Type Modified nucleotides

Formula $C_{50}H_{53}N_9O_{21}P_2$ (free acid)

MW 1178.0 (free acid)

Purity > 95% HPLC

Quantity 0.1 μ mol \approx 0.12mg

Solubility Soluble in water (>3.0mM).

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