

ara-2'-F-NAD⁺

ara-2'-F-NAD⁺ is an analogue of β -NAD⁺ in which the nicotinamide-connected ribose is replaced by 2'-fluoro arabinose. Hydrolysis-resistant slow binding inhibitor of ADP-ribosyl cyclase (i.e. CD 38).

BLG-D148-005 (5 x 0.1 μ mol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.

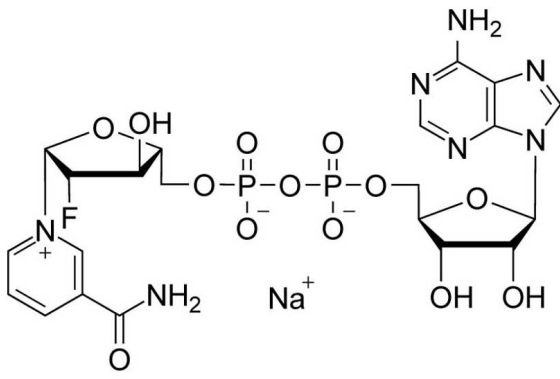
Ordering Information

[Order Online »](#)

BLG-D148-001	0.1 μ mol
BLG-D148-005	5x0.1 μ mol

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	ara-2'-F-NAD ⁺ aliquots (freeze-dried or in solution) are sufficiently stable to be stored up to 7 days at room temperature. Nevertheless, we recommend that the compound should be stored in the freezer, for longer storage periods preferably in freeze-dried form (-20 degrees Celsius necessary, -80 degrees Celsius recommended).
Long Term Storage	-80°C
Shipping	Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	β-ara-2'-Deoxy-2'-fluoro-nicotinamide adenine dinucleotide
Couple Type	Modified nucleotides
Formula	C ₂₁ H ₂₆ FN ₇ O ₁₃ P ₂ (free acid)
MW	665.4 (free acid)
Purity	≥97% (HPLC)
Quantity	0.1 μmol ≈ 67 μg
Solubility	Very well soluble in water and in aqueous solvents (>30mM).

Technical Info / Product Notes λ_{max} = 260nm (ε 20,000, pH 7).



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-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