

2'-Fluoro-NAD⁺

N-2'-F-dAD⁺ is a fluorinated version of the coenzyme nicotinamide adenine dinucleotide (β -NAD⁺). It is a weak inhibitor of human NAD kinase.

*BLG-N073-05 (5 x 1 μ mol pack size) is not sold in the U.S. or Canada.
Please [contact us](#) for available options.*

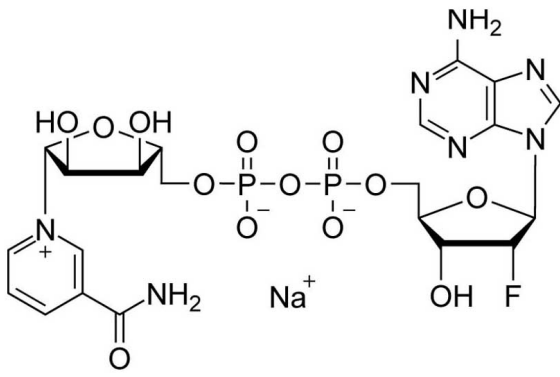
Ordering Information

[Order Online »](#)

BLG-N073-01	1 μ mol
BLG-N073-05	5x1 μ mol

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name N-2'-F-dAD⁺, β-Nicotinamide-2'-deoxy-2'-fluoroadenine dinucleotide . sodium salt

CAS 1514900-83-4

Formula C₂₁H₂₆FN₇O₁₃P₂ (free acid)

MW 665.4 (free acid)

Purity > 95% HPLC



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