

2'-deoxy-NAD⁺

N-2'-dAD is a 2'-deoxyadenosine version of the coenzyme nicotinamide adenine dinucleotide (β -NAD⁺). It is considered to be a substrate for type III CD38 in the enzymatic production of the Transient Receptor Potential Melastatin 2 (TRPM2) channel superagonist dADPR (2'-deoxy-ADPR, Prod. No. BLG-D227). Both N-2'-dAD and dADPR were proved to be present in Jurkat T cells.

BLG-N065-05 (5 x 1 μ mol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.

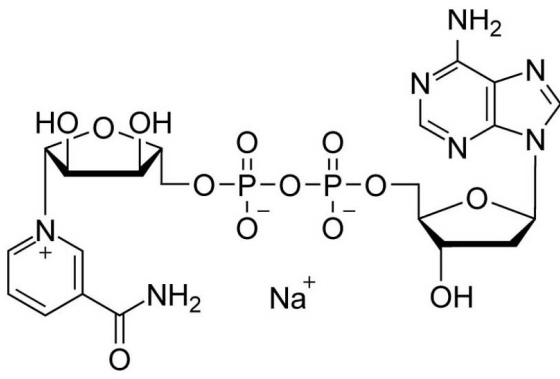
Ordering Information

[Order Online »](#)

BLG-N065-01	1 μ mol
BLG-N065-05	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'-dAD+, β -Nicotinamide-2'-deoxyadenine dinucleotide . sodium salt

CAS 1514900-83-4

Formula $C_{21}H_{27}N_7O_{13}P_2$ (free acid)

MW 674.4 (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