

# ara-2'-F-NAD<sup>+</sup>

ara-2'-F-NAD<sup>+</sup> is an analogue of  $\beta$ -NAD<sup>+</sup> 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  $\mu$ mol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.

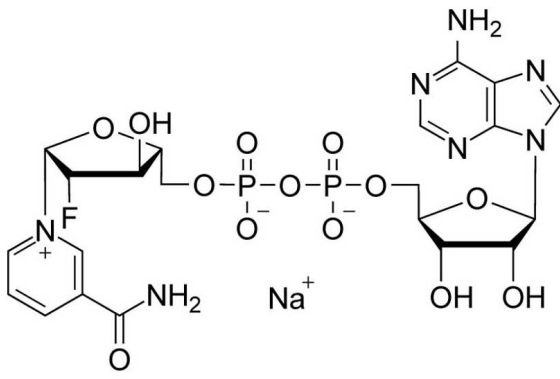
## Ordering Information

[Order Online »](#)

BLG-D148-001	0.1 $\mu$ mol
BLG-D148-005	5x0.1 $\mu$ mol

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

<b>Use/Stability</b>	ara-2'-F-NAD <sup>+</sup> 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).
<b>Long Term Storage</b>	-80°C
<b>Shipping</b>	Ambient Temperature

**Regulatory Status** RUO - Research Use Only

## Product Details

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

**Technical Info / Product Notes** λ<sub>max</sub> = 260nm (ε 20,000, pH 7).



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)