

NAD-Hexylamine, Lithium Salt

A functionalized version of NAD containing an amino group that permits it to be fixable in a gel matrix. This property makes NAD-hexylamine useful for constructing wearable biosensors that comprise a dehydrogenase enzyme such as diaphorase.



Product attributes

CAS number	102029-94-7
------------	-------------

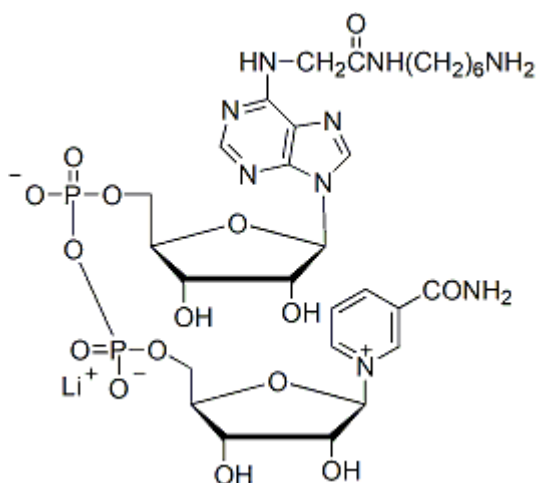
Product Description

NAD (nicotinamide adenine dinucleotide) is a ubiquitous electron carrier and cofactor for a number of oxidoreductase or dehydrogenase enzymes that are central to cellular metabolism.

Biotium's NAD-Hexylamine (N6-((6-aminohexyl)carbamoylmethyl)nicotinamide adenine dinucleotide, or NAD-NH2) is a functionalized version of NAD containing an amino group that permits it to be fixable in a gel matrix. This property makes NAD-hexylamine useful for constructing wearable biosensors that comprise a dehydrogenase enzyme such as diaphorase (1). NAD-Hexylamine may also be used for affinity chromatography, or as a functioning coenzyme to study enzyme mechanisms or topography (2).

NAD exists as a coenzyme redox pair with an oxidized NAD⁺ form and a reduced NADH form. In addition to redox reactions, NAD also participates in ADP-ribose transfer reactions as a donor of ADP-ribose units for enzymes such as ADP-ribosyltransferases and ADP-ribosyl cyclases.

- White solid soluble in water
- Stored desiccated at -20 °C
- C₂₉H₄₂N₉O₁₅Li P₂
- MW: 826
- [102029-94-7]



References

1. JDST(2021) 15, 4, 768-774, [DOI: 10.1177/19322968211008185](https://doi.org/10.1177/19322968211008185)
2. Eur J Bioche (1973) 40, 187, [DOI: 10.1111/j.1432-1033.1973.tb03184.x](https://doi.org/10.1111/j.1432-1033.1973.tb03184.x)

This datasheet was generated on June 4, 2026 at 08:01:03 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/n6-6-aminohexylcarbamoylmethylnicotinamide-adenine-dinucleotide-lithium-salt/>