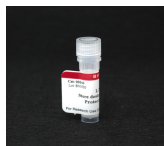


## L-NIO

A non-selective, potent, and irreversible inhibitor of nitric oxide synthase isoforms.



### Product attributes

CAS number	36889-13-1
Molecular weight of antigen	246.1
Storage Conditions	Store at -10 to -35 °C, Desiccate

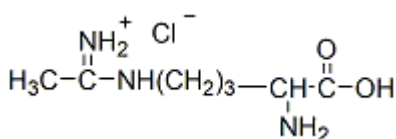
## Product Description

L-NIO (*N*<sup>5</sup>-(1-Iminoethyl)-L-ornithine, dihydrochloride) is a potent, irreversible inhibitor of nitric oxide synthase isoforms.

- Non-selective inhibition of nitric oxide synthase
- White solid soluble in water or MeOH

If a variety of NOS inhibitors are needed, we offer a [NOS Inhibitor Kit](#) which contains four NOS inhibitors: [7-NI \(00240\)](#) (50 mg), [L-NMMA \(00241\)](#) (25 mg), [L-NIL \(00242\)](#) (5 mg) and [L-NIO \(00243\)](#) (5 mg). We also have a selection of dyes and probes for detecting [nitric oxide or reactive oxygen in cells](#).

Molecular structure:



## References

1. Br J Pharmacol, 102(1), 234 (1991), [DOI: 10.1111/j.1476-5381.1991.tb12159.x](https://doi.org/10.1111/j.1476-5381.1991.tb12159.x)
2. Br J Pharmacol, 107(4), 1159 (1992), [DOI: 10.1111/j.1476-5381.1992.tb13423.x](https://doi.org/10.1111/j.1476-5381.1992.tb13423.x)
3. Alcohol Clin Exp Res 33, 1158 (2009), [DOI: 10.1530-0277-2009.00939.x](https://doi.org/10.1530-0277-2009.00939.x)
4. Mol Pain 6, 13 (2010), [DOI: 10.1186/1744-8069-6-13](https://doi.org/10.1186/1744-8069-6-13)
5. Alcohol, 46, 727 (2012), [DOI: 10.1016/j.alcohol.2012.09.001](https://doi.org/10.1016/j.alcohol.2012.09.001)
6. Alcohol, 47, 339 (2013), [DOI: 10.1016/j.alcohol.2013.03.004](https://doi.org/10.1016/j.alcohol.2013.03.004)
7. PLoS ONE, 8(7): e64837. (2013) [DOI: 10.1371/journal.pone.0064837](https://doi.org/10.1371/journal.pone.0064837)
8. Clin Exp Pharmacol Physiol, 41, 246 (2014), [DOI: 10.1111/1440-1681.12207](https://doi.org/10.1111/1440-1681.12207)

This datasheet was generated on June 7, 2026 at 12:55:14 AM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/l-nio-n5-1-iminoethyl-l-ornithine-dihydrochloride/>