

L-NIL

A potent and moderately selective inhibitor of inducible nitric oxide synthase (iNOS) over other NOS isoforms.



Product attributes

CAS number	150403-89-7
Molecular weight of antigen	223.7
Storage Conditions	Store at 2 to 8 °C, Desiccate, Protect from air/store under inert gas

Product Description

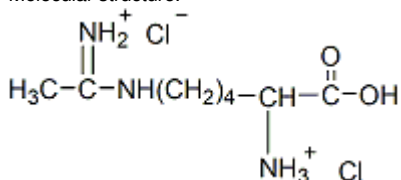
L-NIL (N⁶-(1-Iminoethyl)-lysine, dihydrochloride) is a potent and moderately selective inhibitor of inducible nitric oxide synthase (iNOS).

- Moderately selective inhibition of iNOS over other NOS isoforms
- White solid soluble in water

L-NIL exhibits about 28-fold greater selectivity (IC₅₀ = 3.3 μM) than for the rat brain constitutive enzyme (IC₅₀ = 92 μM). It exhibits about 10-fold greater inhibitory potency than NG-monomethyl-L-arginine in inhibiting γ-interferon-induced NO₂⁻ production (IC₅₀ = 460 nM).

If a variety of NOS inhibitors are needed, we offer a [NOS Inhibitor Kit](#) which contains four 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. J Med Chem 37, 3886 (1994). DOI: [10.1021/jm00049a007](https://doi.org/10.1021/jm00049a007)
2. Eur J Pharmacol 294(2-3), 703 (1995). DOI: [10.1016/0014-2999\(95\)00618-4](https://doi.org/10.1016/0014-2999(95)00618-4)
3. Biochem Pharmacol 51(4), 383 (1996), DOI: [0.1016/0006-2952\(95\)02099-3](https://doi.org/10.1016/0006-2952(95)02099-3)
4. PLoS ONE 8(7): e64837. (2013) DOI: [10.1371/journal.pone.0064837](https://doi.org/10.1371/journal.pone.0064837)

This datasheet was generated on June 10, 2026 at 07:51:14 PM. Visit product page to check for updated information before use.
 Product link: <https://biotium.com/product/l-niin6-1-iminoethyl-lysine-hydrochloride/>