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	223.7
Storage Conditions	Store at 2 to 8 °C, Desiccate, Protect from air/store under inert

Call us: 800-304-5357 Email: btinfo@biotium.com

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 ($IC_{50} = 3.3 \text{ uM}$) than for the rat brain constitutive enzyme (IC₅₀ = 92 uM). 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.

References

- 1. J Med Chem 37, 3886 (1994). <u>DOI: 10.1021/jm00049a007</u>
 2. Eur J Pharmacol 294(2-3), 703 (1995). <u>DOI: 10.1016/0014-2999(95)00618-4</u>
 3. Biochem Pharmacol 51(4), 383 (1996), <u>DOI: 0.1016/0006-2952(95)02099-3</u>
- 4. PLoS ONE 8(7): e64837. (2013) DOI: 10.1371/journal.pone.0064837

This datasheet was generated on November 10, 2025 at 06:42:35 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/l-niln6-1-iminoethyl-lysine-hydrochloride/