



Glowing products for science

Call us : [800-304-5357](tel:800-304-5357)

## 7-Nitroindazole (7-NI)

A selective inhibitor for nitric oxide synthase (NOS) in brain tissues.



### Product attributes

CAS number	2942-42-9
Molecular weight	163.1
Storage Conditions	Store at 2 to 8 °C, Desiccate

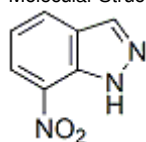
## Product Description

7-Nitroindazole (7-NI) is a selective inhibitor for the brain nitric oxide synthase (NOS) and is often used to study of nitric oxide (NO) activity in the brain.

- Selective inhibition of nitric oxide synthase in brain tissues
- Yellow orange solid soluble in DMSO

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. Meth in Enzymol 268, 393 (1995), [DOI: 10.1016/S0076-6879\(96\)68041-0](#)
2. Stroke 26, 2172 (1995), [DOI: 10.1161/01.STR.26.11.2172](#)
3. Br J Pharmacol 110, 225 (1993), [DOI: 10.1111/j.1476-5381.1993.tb13796.x](#)

This datasheet was generated on January 11, 2026 at 12:07:18 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/7-nitroindazole-7-ni/>