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	

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

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 November 12, 2025 at 09:13:50 AM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/7-nitroindazole-7-ni/