

## Nitric Oxide Generation Kit

This item is discontinued.



### Product attributes

<b>Storage Conditions</b>	See individual components for storage temperature, Desiccate
<b>CAS number</b>	56329-27-2, 136587-13-8, 16142-27-1, 57564-91-7, 79032-48-7

## Product Description

This item is discontinued.

## References

1. J Pharmacol Exp Ther 248, 762 (1989), [J Pharmacol Exp Ther 248, 762 \(1989\)](#)
2. J. Med. Chem 34, 3242 (1991), [DOI: 10.1021/jm00115a013](#)
3. Br J Pharmacol 107, 745 (1992), [DOI: 10.1111/j.1476-5381.1992.tb14517.x](#)
4. Cancer Res. 53, 564-567 (1993), [PMID: 8425188](#)
5. Nature 364, 626 (1993), [DOI: 10.1038/364626a0](#)
6. PNAS 90, 9813 (1993), [DOI: 10.1073/pnas.90.21.9813](#)
7. Neuropharmacology 33, 915 (1994), [DOI: 10.1016/0028-3908\(94\)90190-2](#)
8. Am J Physiol, L9 (1994), [DOI: 10.1152/ajplung.1994.267.1.L9](#)
9. Life Sci 54, 1449 (1994), [DOI: 10.1016/0024-3205\(94\)00600-8](#)
10. Neuron 12, 1235 (1994), [DOI: 10.1016/0896-6273\(94\)90440-5](#)
11. Biochemistry 34, 7177 (1995), [DOI: 10.1021/bi00021a032](#)
12. Biochem Biophys Res Comm 221, 163 (1996), [DOI: 10.1006/bbrc.1996.0563](#)
13. Meth Enzymol 268, 281 (1996), [DOI: 10.1016/S0076-6879\(96\)68030-6](#)
14. Chem. Res. Toxicol. 23, 2 (2010), [DOI: 10.1021/tx900387k](#)
15. Bioorg Med Chem Lett. 21, 6515 (2011), [DOI: 10.1016/j.bmcl.2011.08.071](#)

This datasheet was generated on June 4, 2026 at 08:50:17 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/nitric-oxide-generation-kit/>