

## DEA-NONOate

This item is discontinued.



### Product attributes

CAS number	56329-27-2
Molecular weight of antigen	206.3
Storage Conditions	Store at -10 to -35 °C, Desiccate, Protect from air/store under inert gas

## Product Description

This item is discontinued.

## References

1. Chem. Res. Toxicol. 23, 2 (2010), [DOI: 10.1021/tx900387k](https://doi.org/10.1021/tx900387k)
2. Meth Enzymol 268, 281 (1996), [DOI: 10.1016/S0076-6879\(96\)68030-6](https://doi.org/10.1016/S0076-6879(96)68030-6)
3. J. Med. Chem 34, 3242 (1991), [DOI: 10.1021/jm00115a013](https://doi.org/10.1021/jm00115a013)

This datasheet was generated on June 15, 2026 at 08:10:33 PM. Visit product page to check for updated information before use.  
Product link: <https://biotium.com/product/dea-nonoate-2-nn-diethylamino-diazenolate-2-oxide-diethylammonium-salt/>