

## **DEA-NONOate**

This item is discontinued.



Call us: 800-304-5357

## **Product attributes**

CAS number	56329-27-2
Molecular weight	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), <u>DOI: 10.1021/tx900387k</u>
  2. Meth Enzymol 268, 281 (1996), <u>DOI: 10.1016/S0076-6879(96)68030-6</u>
  3. J. Med. Chem 34, 3242 (1991), <u>DOI: 10.1021/jm00115a013</u>

This datasheet was generated on November 27, 2025 at 09:01:56 PM. Visit product page to check for updated information before use. Product link: <a href="https://biotium.com/product/dea-nonoate-2-nn-diethylamino-diazenolate-2-oxide-diethylammonium-salt/">https://biotium.com/product/dea-nonoate-2-nn-diethylamino-diazenolate-2-oxide-diethylammonium-salt/</a>