Product Information



DPTA NONOate

Item No. 82110

CAS Registry No.: 146724-95-0

Formal Name: (Z)-1-[N-(3-aminopropyl)-N-(3-

ammoniopropyl)amino]diazen-1-ium-1,2-

Synonym: Dipropylenetriamine NONOate

MF: $C_6H_{17}N_5O_2$ FW: 191.2 **Purity:** ≥98%

Stability: ≥1 year at -80°C Supplied as: A crystalline solid λ_{max} : 252 nm UV/Vis.:

Laboratory Procedures

For long term storage, keep dipropylenetriamine NONOate (DPTA NONOate) sealed under nitrogen at -80°C. It should be stable for at least one year. The crystals are sensitive to moisture and become discolored on exposure to air. Keep the vial sealed until use unless your laboratory is equipped with a glove box with an inert atmosphere for the handling of air sensitive compounds.

DPTA NONOate dissociates to the free amine and nitric oxide in a pH-dependent manner following first order kinetics. Alkaline solutions of NONOates (in 0.01 M NaOH) are very stable and can be stored at 0°C for 24 hours. DPTA NONOate is highly soluble in water and relatively concentrated solutions can be prepared for further dilution. To initiate the release of nitric oxide, add a portion of the stock alkaline solution of DPTA NONOate to excess buffer of pH 7.0-7.4. The half-life of DPTA NONOate is 5 hours at 22°C in 0.1 M phosphate buffer (pH 7.4). The decomposition of NONOates is nearly instantaneous at pH 5.1

DPTA NONOate is a NO donor. It spontaneously dissociates in a pH-dependent, first-order process with a half-life of three hours and five hours at 37°C and 22-25°C, pH 7.4, respectively, to liberate 2 moles of NO per mole of parent compound.1,2

References

- Hrabie, J.A., Klose, J.R., Wink, D.A., et al. New nitric oxide-releasing zwitterions derived from polyamines. J. Org. Chem. 58, 1472-1476 (1993).
- Keefer, L.K., Nims, R.W., Davies, K.M., et al. 'NONOates' (1-substituted diazen-1-ium-1,2-diolates) as nitric oxide donors: Convenient nitric oxide dosage forms. Methods Enzymol. 268, 281-293 (1996).

WARNING: This product is for laboratory research only: not for administration to humans. Not for human or veterinary DIAGNOSTIC OR THERAPEUTIC USE.

MATERIAL SAFETY DATA

MALERIAL SAFELY DALA
This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. Before use, the user must review the complete Material Safety Data Sheet, which has been sent via email to your institution

WARRANTY AND LIMITATION OF REMEDY

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular bose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman warrants only to the original customer that the material will meet our specifications

Layman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have any obligation or liability, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees.

Buyer's exclusive remedy and Cayman's sole liability hereunder shall be limited to a refund of the purchase price, or at Cayman's option, the replacement, at no cost to Buyer, of all material that

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within rty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our Warranty and Limitation of Remedy located on our website and in our catalog.

Copyright Cayman Chemical Company, 06/19/2012

Phone

(734) 971-3640

(800) 364-9897 (734) 971-3335

custserv@caymanchem.com

www.caymanchem.com

Cayman Chemical

Mailing address 1180 E. Ellsworth Road Ann Arbor, MI 48108 USA