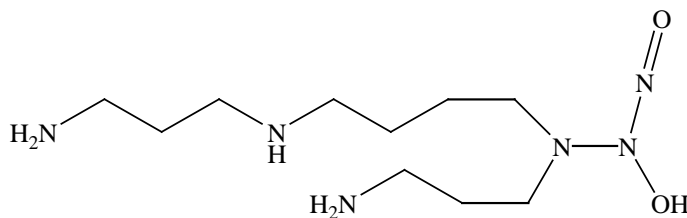


## PRODUCT AND SAFETY DATA SHEET

**PRODUCT NAME:** Spermine NONOate**CATALOG #:** 00224**MOLECULAR INFORMATION:** C<sub>10</sub>H<sub>26</sub>N<sub>6</sub>O<sub>2</sub>  
Mwt: 262.4  
[136587-13-8]**PROPERTIES:**

<b>Color &amp; Form</b>	Crystalline solid
<b>Purity</b>	≥ 98%
<b>Solubility</b>	Soluble in water
<b>Absorption</b>	252nm
<b>Extinction Coefficient</b>	8,500

**STORAGE AND HANDLING:**

Hygroscopic. Store desiccated at -20 °C. For long term storage, keep sealed at -80 °C, and it will be stable for at least one year

**APPLICATION:**Spermine NONOate dissociates to the free amine and nitric oxide in a pH-dependent manner following first order kinetics. To initiate the release of nitric oxide, add a portion of the stock alkaline solution of Spermine NONOate to excess buffer of pH 7.0-7.4. The half-life of Spermine NONOate is 39 minutes at 37°C in 0.1 M phosphate buffer (pH 7.4). The decomposition of the NONOate is nearly instantaneous at pH 5.<sup>1</sup>Ref: 1) *J. Med. Chem* **34**, 3242-3247 (1991); 2) *Cancer Res.* **53**, 564-567 (1993); 3) *Proc. Natl. Acad. Sci. USA* **90**, 9813-9817 (1993)**TOXICITY:** Unknown

<b>FIRST AID:</b>	Potentially harmful; avoid prolonged or repeated exposure. Avoid getting in eyes, on skin, or on clothing. Wash thoroughly after handling. If eye or skin contact occurs, wash affected areas with plenty of water for 15 minutes and seek medical advice. In case of inhaling or swallowing, move individual to fresh air and seek medical advice immediately.
-------------------	---

<b>Disclaimer:</b> Materials from Biotium are sold for research use only, and are not intended for food, drug, household, or cosmetic use. Biotium is not liable for any damage resulting from handling or contact with this product.
---