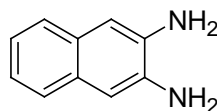


PRODUCT AND SAFETY DATA SHEET

PRODUCT NAME: 2,3-Diaminonaphthalene**CATALOG #:** 00301**MOLECULAR INFORMATION:** C₁₀H₁₀N₂
Mwt: 158**PROPERTIES:**

Color & Form	Brown Crystalline solid
Purity	≥ 98% by HPLC
Solubility	Soluble in DMSO
Absorption	N/A
Extinction Coefficient	N/A

STORAGE AND HANDLING: Store desiccated at 4°C**APPLICATION:** 2,3-Diaminonaphthalene reacts with nitrosonium, which is formed from NO, to form the fluorescent dye 1 *H*-naphthotriazole.^{1,2} This method can be used to detect 10 nM to 10 μM of nitrite (NO₂⁻) and is compatible with 96-well format.³Ref.: 1) *Luminescence* **14**, 283 (1999); 2) *Methods Enzymol.* **268**, 105(1996); 3) *Anal. Biochem.* 214, 11(1993)**TOXICITY:** Unknown

FIRST AID:	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.
-------------------	---

Disclaimer: <i>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.</i>
--