

**2,3-Diaminonaphthalene \*CAS 771-97-1\***Catalog number: 15221  
Unit size: 100 mg**Product Details**

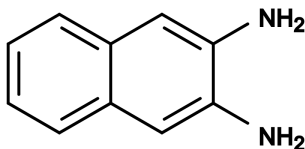
---

Storage Conditions	Freeze (<-15 °C), Minimize light exposure
Expiration Date	12 months upon receiving

**Chemical Properties**

---

Appearance	Light brown solid
Molecular Weight	158.20
Soluble In	DMSO
Chemical Structure	

**Applications**

---

The role of nitric oxide (NO) as a putative mediator of neuronal death can be understood best if NO is detected directly. 1,2-Diaminonaphthalene (DAN) is a fluorogenic probe that is used to detect nitric oxide (NO) productions in live cell and animals. The reaction product of DAN with NO is a triazole with blue fluorescence. The DAN-based methodology of NO assessment is convenient and applicable to numerous living systems both in vivo and in vitro. As low as 10 nM nitrite can be detected using DAN.