

**NBD-F [4-Fluoro-7-nitrobenzofurazan] \*CAS 29270-56-2\***

Catalog Number: 821

Unit Size: 5 mg

**Product Details**

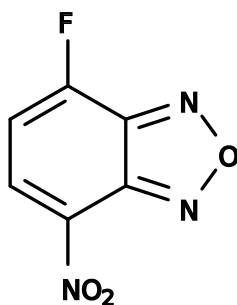
---

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

**Chemical Properties**

---

Appearance	Solid
Molecular Weight	183.1
Soluble In	DMSO
Chemical Structure	

**Spectral Properties**

---

Excitation Wavelength	467 nm
Emission Wavelength	539 nm

**Applications**

---

NBD-F has properties and applications similar to those of NBD-Cl. Upon reacting with amines and thiol compounds it generates the same fluorescent adducts as those of NBD-Cl. Compared with NBD-Cl, NBD-F is much more reactive, and should be more carefully stored. For example, the reaction of NBD fluoride with glycine is reported to be 500 times faster than the reaction of NBD chloride with glycine. Both NBD chloride and NBD fluoride are extensively used as derivatization reagents for chromatographic analysis of amino acids and other low molecular weight amines.