## **NBD Fluoride**

NBD fluoride (4-Fluoro-7-nitrobenzo-2-oxa-1,3-diazole) is similar to NBD chloride (90045) except that the former is much more reactive



### **Product attributes**

Excitation/Emission

465/535 nm (reaction product with primary amine); 485/540 nm (reaction product with secondary amine)

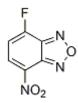
Call us: 800-304-5357 Email: btinfo@biotium.com

# toward amines.

## **Product Description**

NBD fluoride (4-Fluoro-7-nitrobenzo-2-oxa-1,3-diazole) is similar to NBD chloride (90045) except that the former is much more reactive toward amines. Both reagents are nonfluorescent until they react with primary or secondary amines to produce a fluorescent product. NBD chloride and NBD fluoride have been extensively used as derivatizing reagents for chromatography analysis of amino acids and low molecular weight amines.

- λEx/λ Em (MeOH) = 465/535 nm (for reaction product with primary amines); 485/540 nm (for reaction product with secondary amines)
- Yellow solid soluble in DMF
- Store at -20°C and protect from light
- C<sub>6</sub>H<sub>2</sub>FN<sub>3</sub>O<sub>3</sub>
- MW:183



## References

- 1. Anal Chim Acta 130, 377 (1981).
- 2. Biochem J 108, 155 (1968).
- 3. Anal Biochem 116, 471 (1981).
- 4. Anal Chim Acta 290, 3 (1994).

This datasheet was generated on November 4, 2025 at 02:05:55 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/nbd-fluoride-4-fluoro-7-nitrobenzo-2-oxa-13-diazole/