

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 toward amines.



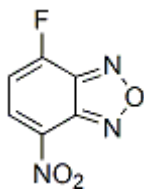
Product attributes

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

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.

- $\lambda_{Ex}/\lambda_{Em}$ (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
- $C_6H_2FN_3O_3$
- 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 June 4, 2026 at 07:56:51 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/>