

NBD Chloride

NBD chloride (4-Chloro-7-nitrobenzo-2-oxa-1,3-diazole) is nonfluorescent until it reacts with primary or secondary amines to produce a fluorescent product. NBD chloride has been extensively used as a derivatizing reagent for chromatography analysis of amino acids and low molecular weight amines.



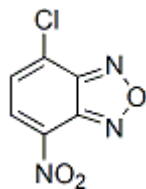
Product attributes

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

Product Description

NBD chloride (4-Chloro-7-nitrobenzo-2-oxa-1,3-diazole) is nonfluorescent until it reacts with primary or secondary amines to produce a fluorescent product. NBD chloride has been extensively used as a derivatizing reagent for chromatography analysis of amino acids and low molecular weight amines.

- $\lambda_{Ex}/\lambda_{Em}(\text{MeOH}) = 465/535 \text{ nm}$ (for reaction product with primary amines); $485/540 \text{ nm}$ (for reaction product with secondary amines)
- Yellow solid soluble in DMF
- Store at -20°C and protect from light
- $\text{C}_6\text{H}_2\text{ClN}_2\text{O}_3$
- MW: 199.55



References

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

This datasheet was generated on December 30, 2025 at 02:51:43 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/nbd-chloride-4-chloro-7-nitrobenzo-2-oxa-13-diazole/>