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.

atictum 680

Product attributes

Excitation/Emission

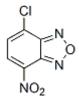
465/535 nm (reaction product with primary amine); 485/540 nm (reaction product with secondary amine)

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

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.

- λ_{Ex}/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
- C₆H₂CIN₃O₃
- 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 November 3, 2025 at 12:40:20 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/