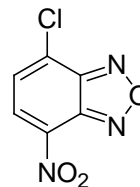


PRODUCT AND SAFETY DATA SHEET

PRODUCT NAME: NBD Chloride (4-chloro-7-nitrobenzo-2-oxa-1,3-diazole)**CATALOG #** 90045**MOLECULAR INFORMATION:** C₆H₂ClN₂O₃
MWt: 199.55**PROPERTIES:**

Color & Form Yellow solid
Purity ≥ 98% by TLC
Solubility Soluble in DMF
Absorption/Emission 465/535 nm (for reaction product with primary amines); 485/540 nm (for reaction product with secondary amines)

Extinction Coefficient

STORAGE AND HANDLING: Store at -20 °C and protect from light

APPLICATION: NBD chloride 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.

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

TOXICITY: Not established. Not listed by NTP, IARC or OSHA.

FIRST AID:	Potentially harmful. Avoid prolonged or repeated exposure. Avoid getting in eyes, on skin, or on clothing. Wash thoroughly after handling. If eye or skin contact occurs, wash affected areas with plenty of water for 15 minutes and seek medical advice. In case of inhaling or swallowing, move individual to fresh air and seek medical advice immediately.
-------------------	---

Disclaimer: <i>Materials from Biotium are sold for research use only, and are not intended for food, drug, household, or cosmetic use. Biotium is not liable for any damage resulting from handling or contact with this product.</i>
--