

## NBD-Ethylenediamine

The fluorescence of NBD is highly dependent on the environment; it is virtually nonfluorescent in water but becomes strongly green fluorescent in hydrophobic media. NBD-ethylenediamine can be readily coupled to activated carboxy groups or sulfonyl chlorides.



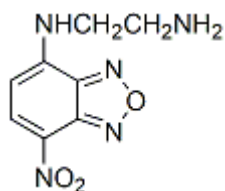
### Product attributes

Excitation/Emission	466/535 nm
---------------------	------------

## Product Description

The fluorescence of NBD is highly dependent on the environment; it is virtually nonfluorescent in water but becomes strongly green fluorescent in hydrophobic media. NBD-ethylenediamine can be readily coupled to activated carboxy groups or sulfonyl chlorides.

- $\lambda_{Ex}/\lambda_{Em}$  (MeOH) = 466/535 nm
- Yellow solid soluble in DMSO or DMF
- Store at 4 °C and protect from light
- $C_8H_9N_5O_3$
- MW: 223.19



This datasheet was generated on December 28, 2025 at 02:51:39 AM. Visit product page to check for updated information before use.  
Product link: <https://biotium.com/product/nbd-ethylenediamine/>