

ABD-F

ABD-F is a useful reagent for the sensitive and specific detection of thiols both by TLC and HPLC (1,2). When combined with a reducing agent such as tributylphosphine, it can also be used for the determination of disulfides in peptides and proteins (3,4).



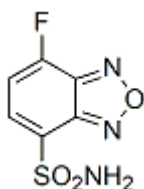
Product attributes

Excitation/Emission	375/508 nm
---------------------	------------

Product Description

ABD-F is a useful reagent for the sensitive and specific detection of thiols both by TLC and HPLC (1,2). When combined with a reducing agent such as tributylphosphine, it can also be used for the determination of disulfides in peptides and proteins (3,4).

- $\lambda_{Ex}/\lambda_{Em}$ (N-acetylcysteine conjugate at pH 2) = 375/508 nm
- Of-white solid soluble in DMSO or DMF
- Store desiccated at 4 °C and protect from light especially in solution
- $C_7H_7FN_3O_3S$
- MW: 232



References

1. Anal Chem 56, 2461 (1984).
2. J Pharm Biomed Anal 24, 1137 (2001).

This datasheet was generated on June 17, 2026 at 05:53:53 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/abd-f/>