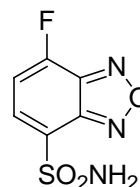


PRODUCT AND SAFETY DATA SHEET

PRODUCT NAME: ABD-F (4-(aminosulfonyl)-7-fluoro-2,1,3-benzoxadiazole)**CATALOG #:** 91046**MOLECULAR INFORMATION:** C₆H₄FN₃O₃S
Mwt: 217.2
[91366-65-3]**PROPERTIES:**

Color & Form	Off-white solid.
Purity	≥ 98% by TLC
Solubility	Soluble in DMSO or DMF
$\lambda_{ex} \backslash \lambda_{em}$	(N-acetylcysteine conjugate at pH 2) = 375/508 nm
Extinction Coefficient	N/A

STORAGE AND HANDLING:

Store desiccated at ≤4 °C and protect from light especially when in solution.

APPLICATION: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}Ref.: 1) *Anal. Chem.* **56**, 2461(1984); 2) *J. Pharm. Biomed. Anal.* **24**, 1137(2001); 3) *Anal. Biochem.* **214**, 128(1993); 4) *J. Biol. Chem.* **264**, 7185(1989)**TOXICITY:**

Unknown.

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: 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.