

# PRODUCT DATA SHEET

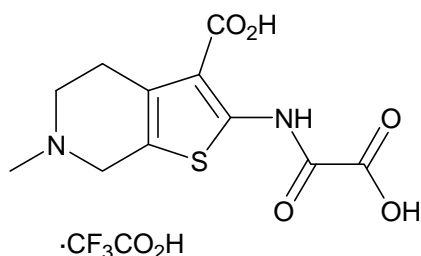


PRODUCT: BML-267  
6-Methyl-2-(oxalylamino)-4,5,6,7-tetrahydrothieno[2,3-c]pyridine-3-carboxylic acid trifluoroacetic acid salt

CATALOG NO.: PR-121

LOT NO.: temp

STRUCTURE:



PHYSICAL APPEARANCE: White solid

MOLECULAR FORMULA:  $C_{11}H_{12}N_2O_5S \cdot C_2HF_3O_2$  MOLECULAR WEIGHT: 398.3

PURITY: 95% (NMR)

SOLUBILITY: Soluble in: DMSO 10 mg/mL

STORAGE: Store as supplied, at  $-20^{\circ}C$  for up to 1 year. Store solutions at  $-20^{\circ}C$  for up to 1 month.

APPLICATION NOTES: Potent and selective protein tyrosine phosphatase 1B (PTP1B) inhibitor ( $IC_{50}=11 \mu M$ )<sup>1</sup>. A similar N-des-methyl analog has been reported<sup>2</sup> and has been used to probe the mechanism of PTP1B-mediated inhibition of leptin signaling<sup>3</sup>. BML-267 and analogs have limited cell permeability<sup>3</sup>. A cell permeable diethyl ester prodrug of this compound is available (catalog # PR-129).

REFERENCES: 1. H.S. Andersen *et al. J. Med. Chem.* 2002 **45** 4443  
2. L.F. Iversen *et al. J. Biol. Chem.* 2000 **275** 10300  
3. I.K. Lund *et al J. Mol. Endocrinol.* 2005 **34** 339

**The pharmacological and toxicological properties of this product have not been fully investigated. Exercise caution in use and handling. This product must not be used in humans.**