

## Data Sheet

### MTAP

human, recombinant, N-terminal GST tag  
 Catalog # 50305  
 Lot#: 80902                      Conc.: 2.5 mg/ml

**Formulated in:** 45 mM Tris-HCl, pH 8.0,  
 124 mM NaCl, 2.4 mM KCl, 18 mM  
 glutathione, 10% glycerol, and 3 mM DTT.

**Stability:** >6 months at -80°C

**References:**

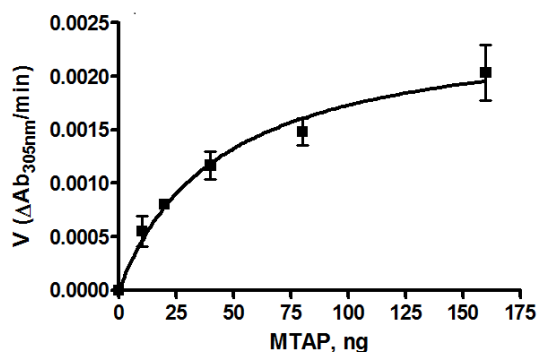
1. Fernandez-Irigoyen, J. *et al.*, *Biochem. J.* **411** (2), 457-465 (2008)
2. Hellerbrand, C. *et al.*, *Carcinogenesis* **27** (1), 64-72 (2006)
3. Ong, S-E., *et al.* *PNAS* **106**(12): 4617-4622 (2009)

**Description:** Human MTAP (GenBank Accession No. NM\_002451), amino acids 2-end, with N-terminal GST tag, MW=57 kDa, expressed in an *E. coli* expression system.

**Assay Conditions:** MTAP activity was measured using a xanthine oxidase-coupled method. MTA cleavage by MTAP produce adenine, which is then converted to 2,8-dihydroxyadenine by xanthine oxidase, leading to an increase in absorbance at 305 nm. 50 uL reaction mixtures contained MTAP, 50 mM KPO<sub>4</sub>, pH 7.4, 200 uM MTA, 0.02 units Xanthine Oxidase, and 1 mM DTT. Activity was measured at 37°C by continuous monitoring the absorbance at 305 nm for 30 min.

**Application:** Suitable for protein binding studies, Western blotting, or as an immunogen.

### Quality Assurance



**10% SDS-PAGE  
 Coomassie staining**

**Lane 1:**  
 4 µg MTAP  
**Lane 2:**  
 Protein Marker

**MW:** 57 kDa  
**Purity:** ≥80%

