

## Data Sheet

### **MAT2A, His-Tag**

Human, Recombinant, N-terminal His-Tag

**Catalog #:** 71401

**Lot #:** 171222-1      **Conc.:** 3.2 mg/ml

**Formulated in:** 40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween, 20% glycerol, additional 250mM imidazole

**Stability:** At least 6 months at  $-80^{\circ}\text{C}$ . Avoid freeze/thaw cycles. Protein may be diluted to  $\geq 100 \mu\text{g/ml}$  in PBS + glycerol and stored at  $-80^{\circ}\text{C}$ .

**References:**

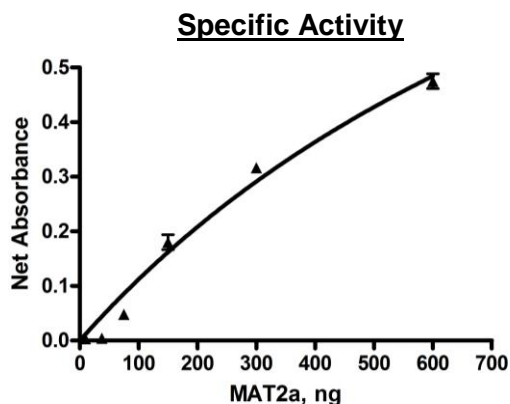
1. Latasa, M., et al. *The FASEB Journal*. 2001; **15(7)**: 1248-1250.
2. Liu, Q., et al., *Molecular Cancer Therapeutics*. 2011; **10(6)**:1113-23.

**Description:** Human MAT2A, also known as methionine adenosyltransferase II alpha, MAT2, and adoMet synthase 2. GenBank Accession No. NM\_005911, a.a. 2-395 (end) with an N-terminal His-tag, expressed in an E. coli cell expression system. MW = 45 kDa.

**Assay Description:** MAT2A #71401 was tested according to the MAT2A Inhibitor Screening Assay Kit catalog #71402 with enzyme tested from 0.9375 ng/ $\mu\text{l}$ -120 ng/ $\mu\text{l}$ . The reactions were conducted at room temperature for 45 minutes in a 25  $\mu\text{l}$  mixture containing 50 mM Tris-HCl, pH 8.0, 10 mM  $\text{MgCl}_2$ , 50 mM KCl, 0.1 mg/ml BSA, various amount of MAT2a enzyme, 200  $\mu\text{M}$  L-Methionine, and 200  $\mu\text{M}$  ATP. After the enzymatic reactions, 50  $\mu\text{l}$  of colorimetric detection reagent was added to each well and the reactions were incubated for another 15 min before the absorbance at 630 nm of each well was measured using a Tecan Infinite M1000 microplate reader.

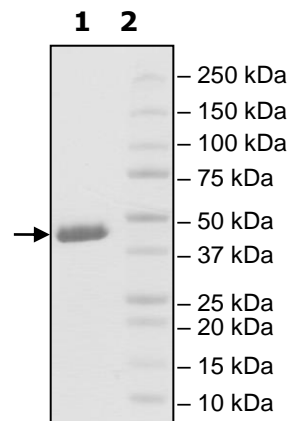
**Application:** Useful for studying protein binding and screening small molecules.

### Quality Assurance



**4-20% SDS-PAGE  
Coomassie staining**

**Lane 1:**  
MAT2A  
**Lane 2:**  
Protein Marker  
**MW:** 45 kDa  
**Purity :**  $\geq 90\%$



OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**  
 Or you can Email us at: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)  
 Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)