

ABAT (4-aminobutyrate aminotransferase)

**CATALOG NO.:** MSC-11-676

**DESCRIPTION:** Human recombinant ABAT (residues 29-500; Genbank Accession # NM\_020686.6; MW = 64.8 kDa) expressed in *E. coli* with an N-terminal His-tag and C-terminal StrepII tag. Catalyzes the neurotransmitter gamma-aminobutyric acid (GABA) into succinic semialdehyde.

**PURITY:** > 95% by SDS-PAGE.

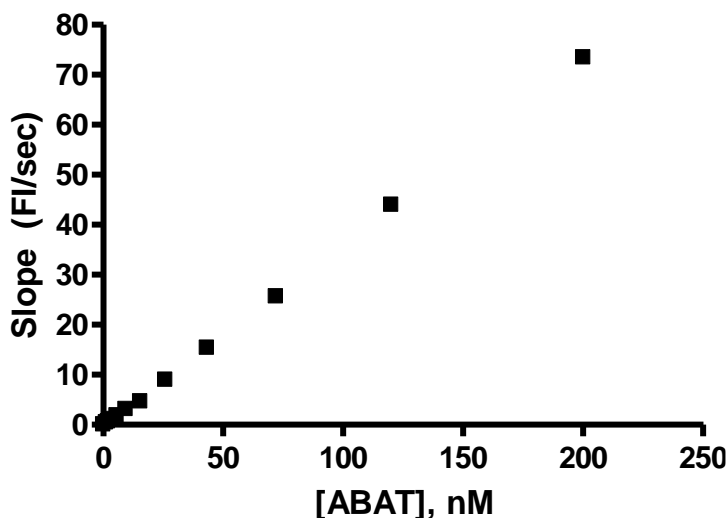
**SUPPLIED AS:** \_\_\_ µg/µL in 50 mM Tris, pH 7.5, 500 mM NaCl, 0.5 mM TCEP, 10% (w/v) glycerol as determined by OD280.

**ASSAY CONDITIONS:** RBC's ABAT displays activity in a succinate dehydrogenase and diaphorase/resazurin coupled assay in the presence of 5 mM GABA and 2 mM aKG, along with detection mixture of 0.2 mM SSDH, 0.5 mM NAD, 0.015 mg/mL diaphorase and 50 µM resazurin. Reaction conditions are: 50 mM potassium pyrophosphate, pH 8.5, 0.01% Triton X-100, substrates at concentrations indicated above.

**STORAGE:** -70°C. Thaw quickly and store on ice before use. The remaining, unused, undiluted protein should be snap frozen, for example in a dry ice/ethanol bath or liquid nitrogen. Minimize freeze/thaws if possible, but very low volume aliquots (<5 µl) or storage of diluted enzyme is not recommended.



Coomassie blue stained SDS-PAGE (4-20% acrylamide) of 4 µg of RBC ABAT. MW markers (left) are, from top, 220, 160, 120, 100, 90, 80, 70, 60, 50, 40, 30, 25, 20, 15, 10 kDa.



**ABAT activity assay.** Activity of RBC's ABAT was monitored using a succinate dehydrogenase/diaphorase/resazurin coupled assay. Variable amounts of ABAT were incubated with 5 mM GABA and 2 mM aKG along with detection mixture of 0.2 µM SSDH, 0.5 mM NAD, 0.015 mg/mL diaphorase and 50 µM resazurin. CLARIOstar plate reader (BMG Labtech) was used for real time monitoring with Ex/Em 530/595nm.

This product is not intended for therapeutic or diagnostic use in animals or in humans.

Reaction Biology

1 Great Valley Parkway, Malvern PA, USA 19355

requests@reactionbiology.com www.reactionbiology.com