

Data Sheet

PSAT1

Human, recombinant, N-terminal GST tag

Catalog #: 71085

Lot #: 130117 Conc.: 2.44 mg/ml

Formulated in: 40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 16 mM glutathione, 0.04% Tween-20, and 20% glycerol.

Stability: >6 months at -80°C

Reference(s):

1. Baek, J.Y., *et al.*, *Biochem J.* 2003 Jul 1;**373(Pt 1)**: 191-200.
2. Hart, C.E., *et al.*, *Am J Hum Genet.* 2007 May;**80(5)**:931-7.

Description:

Human phosphoserine aminotransferase 1 (PSAT1), also known as EPIP, amino-acids 2 – 370 (end) (GenBank Accession No. NM_058179) with N-terminal GST tag, MW = 67.1 kDa, expressed in an *E. coli* expression system.

Application:

Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

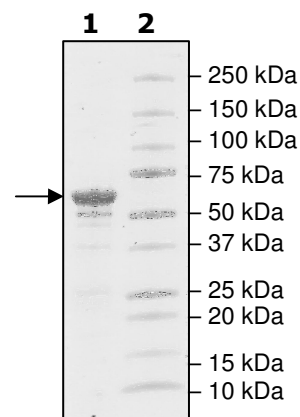
Quality Assurance

4-20% SDS-PAGE Coomassie staining

Lane 1:
3.6 µg PSAT1

Lane 2:
Protein Marker

MW: 67.1 kDa
Purity: ≥76%



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

Please visit our website at: www.bpsbioscience.com



6044 Cornerstone Court West, Suite E
San Diego, CA 92121
Tel: 1.858.829.3082
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Sequence:

MSPILGYWKIKGLVQPTRLLEYLEEKYEEHLYERDEGDKWRNKKFELGLEFPNLPYYI
DGDVKLTQSMARIYIADKHNMLGGCPKERAIEISMLEGAVLDIRYGVSRAYSKDFETLK
VDFLSKLPEMLKMFEDRLCHKTYLNGDHVTHPDFMLYDALDVVLYMDPMCLDAFPKLV
CFKKRIEAIQIDKYLKSSKYIAWPLQGQWQATFGGGDHPPKSDLEVLFQGPLGSM
DAP
RQVVNFGPGPAKLPHSVLLEIQKELLDYKGVGISVLEMSHRSSDFAKIINNTENLVREL
LAVPDNYKVIFLQGGGCGQFSVPLNLIGLKAGRCADYVVTGAWSAKAAEEAKKFGT
INIVHPKLGSYTKIPDPSTWNLNPDASYVYYCANETVHGVEFDVIPDVKGAVLVCDMSS
NFLSKPVDVSKFGVIFAGAQNVSAGVTVVIVRDDLLGFALRECPSVLEYKVQAGNS
SLYNTPPCFSIYVMGLVLEWIKNNGGAAAMEKLSSIKSQTIEIIDNSQGFYVCPVEPQ
NRSKMNIPFRIGNAKGDDALEKRFLDKALELNMLSLKGHRVGGIRASLYNAV TIEDV
QKLA AFMKKFLEMHQL

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
Please visit our website at: www.bpsbioscience.com