

## Data Sheet

### **ATAT1 (2 – 224)**

Human, recombinant, N-terminal GST-tag  
Catalog #: 50072  
Lot#: 120410-RT      Conc.: 0.14 mg/ml

**Formulated in:** 40 mM Tris-HCl, pH 8.0,  
110 mM NaCl, 2.2 mM KCl, 16 mM  
Glutathione, 20% glycerol, and 3 mM DTT.

**Stability:** >6 months at –80 °C

**References:**

1. Akella, J.S., *et al.*, *Nature*. 2010 Sep 9;  
**467(7312)**: 218 – 222.
2. Lloyd, R.A., *et al.*, *Anal. Biochem.* 1994  
Jan ; **216(1)** : 42 -46.

**Description:**

Human alpha tubulin acetyltransferase 1,  
or ATAT1, GenBank Accession No.  
NM\_001031722 (var1), a.a. 2 – 224, with  
N-terminal GST-tag, MW=52 kDa,  
expressed in an *E. coli* expression  
system.

**Application:**

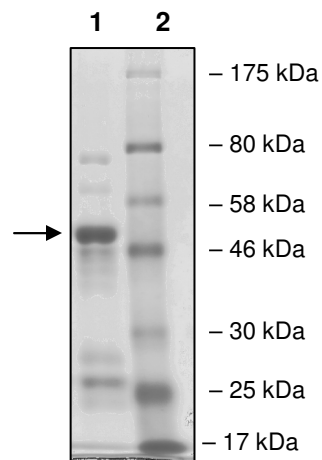
Suitable for protein binding studies as an  
immunogen or as a protein standard.

## ***Quality Assurance***

**10% SDS-PAGE**  
**Coomassie staining**

**Lane 1:**  
2.1 µg ATAT1  
**Lane 2:**  
Protein Marker

**MW:** 52 kDa  
**Purity:** ≥40%



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)



6044 Cornerstone Court W, Ste E  
San Diego, CA 92121  
**Tel:** 1.858.829.3082  
**Fax:** 1.858.481.8694  
**Email:** [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

**Protein Sequence:**

MSPILGYWKIKGLVQPTRLLEYLEEKYEEHLYERDEGDKWRNKKFELGLEFPNLPYYI  
DGDVKLTQSMARIYIADKHNMLGGCPKERAISMLEGAVLDIRYGVSRAYSDFETLK  
VDFLSKLPEMLKMFEDRLCHKTYLNGDHVTHPDFMLYDALDVVLYMDPMCLDAFPKLV  
CFKKRIEAIQIDKYLKSSKYIAWPLQGWQATFGGGDHPPKSDLEVLFGGPLGSWLTW  
PFCFLTITLREEGVCHLESVDLQQQIMTIIDELGKASAKAQNLSAPITSASRMQSNRHV  
VYILKDSSARPAGKGAIGFIKVGKLFVLDDREAHNEVEPLCILDFYIHESVQRHGHG  
RELQYMLQKERVEPHQLAIDRPSQKLLKFLNKHYNLETTVPQVNNFVIFEGFFAHQH  
RPPAPSLRATRHSRAAAVDPTPAAPARKLPPKRAEGDIKPYS

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)