

Data Sheet

HSP90 α , His-Avi-Tag, Biotin-labeled (Mouse)

Mouse, recombinant, N-terminal His-Avi-tag,
Biotin-labeled

Catalog #: 100751

Lot#: 200826 **Conc.:** 3.44 mg/ml

Formulated in: 40 mM Tris-HCl, pH 8.0,
110 mM NaCl, 2.2 mM KCl, 0.04%
Tween20, 20% glycerol, 3 mM DTT.

Stability: At Least 6 months at -80°C .
Avoid freeze/thaw cycles.

References:

- Wayne N., *et al.* *J Biol Chem.* 2007 Nov 30;**282(48)**:35386-95.
- Pearl L.H., *et al.* *Biochem J.* 2008 Mar 15;**410(3)**:439-53.

Description: Mouse Heat Shock Protein 90 alpha (HSP90 α), GenBank Accession No. NM_010480.5, a.a. 535-733(end) with N-terminal His-Avi-tag, expressed in an *E. coli* expression system, MW=25 kDa. This protein runs at a higher by SDS-PAGE due to glycosylation. This protein is enzymatically biotinylated using Avi-Tag™ technology. Biotinylation is confirmed to be $\geq 90\%$

Quality Assurance

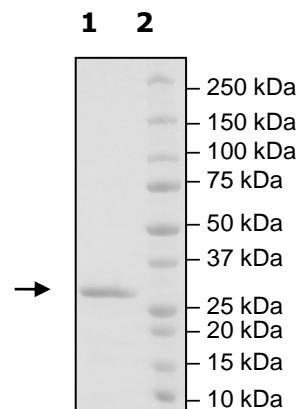
**4-20% SDS-PAGE
Coomassie staining**

Lane 1:
HSP90a

Lane 2:
Protein Marker

Purity: >90%

MW: 25 kDa



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

To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**
Or you can Email us at: info@bpsbioscience.com
Please visit our website at: www.bpsbioscience.com