

Data Sheet

HSP90 β (C-terminal), Biotin-labeled

Human, recombinant, N-terminal His-tag,
N-terminal Avi-tag™

Catalog #: 50313

Lot#: 130607-G **Conc.:** 0.68 mg/ml

Formulated in: 40 mM Tris-HCl pH 8, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween-20, 20% glycerol

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

References:

1. Milicevic, Z. *et al.*, *Int. J. Oncol.* 2008 Jun;**32(6)**:1169-78.
2. Chao, C.C. *et al.*, *J. Cell Biochem.* 2008 Jul 1;**104(4)**:1286-96.

Description:

Human Heat Shock Protein 90 beta (HSP90 β), GenBank Accession No. NM_001271969, a.a. 527-724 (end) with N-terminal His-tag and N-terminal Avi-tag™, MW = 25 kDa, expressed in an *E. coli* expression system and enzymatically biotin-labeled using Avi-tag™ technology.

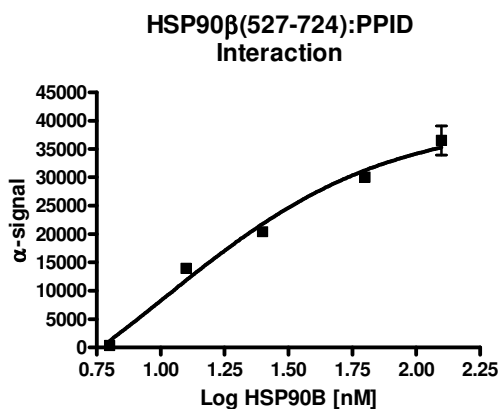
Assay Conditions:

HSP90 β (C-terminal):PPID binding activity measured using HSP90 β (C-terminal) Inhibitor Screening Assay Kit, BPS Bioscience, Catalog # 50314.

Application:

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

Quality Assurance

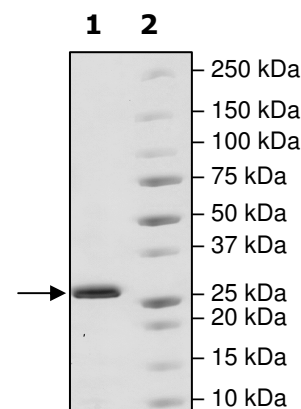


4-20% SDS-PAGE Coomassie staining

Lane 1:
2 μg HSP90 β

Lane 2:
Protein Marker

MW: 25 kDa
Purity: $\geq 90\%$



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

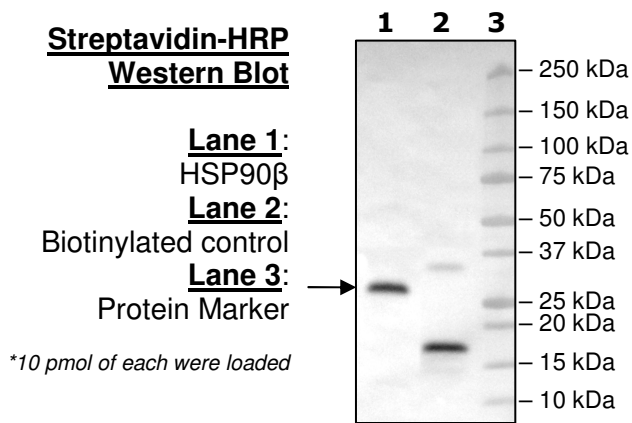
To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**

Or you can Email us at: info@bpsbioscience.com

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:

MHHHHHGGGLNDIFEAQKIEWHEEFDGKSLVSVTKEGLELPEDEEEKKKMEESKAK
FENLCKLMKEILDKKVEKVTISNRLVSSPCCIVTSTYGWTANMERIMKAQALRDNSTM
GYMMAKKHLEINPDHPIVETLRQKAEADKNDKAVKDLVLLFETALLSSGFSLEDPQT
HSNRIYRMIKLGLGIDEDEVAAEENAAVPDEIPPLEGDEDASRMEEVD

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