

Data Sheet

Aha1, His-Tag

Human, recombinant, C-terminal His-Tag

Catalog #: 50291

Lot#: 130430-250E1 **Conc.:** 1.24 mg/ml

Formulated in: 40mM Tris pH8, 110mM NaCl, 2.2mM KCl, 3mM DTT, 20% glycerol, 200mM Imidazole.

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

References:

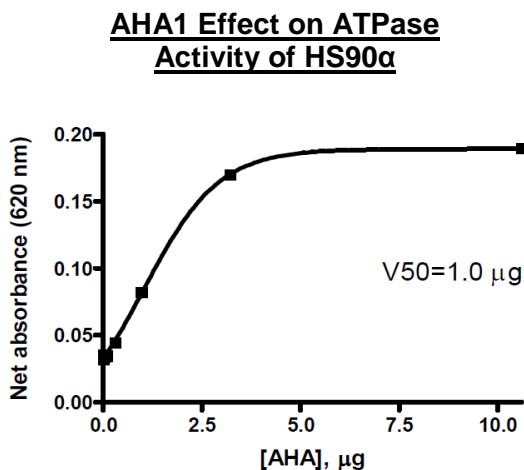
1. Wang, X. *et al.*, *Cell* **127** (4), 803-815 (2006).
2. Lotz, G.P. *et al.*, *J. Biol. Chem.* **278** (19), 17228-17235 (2003).

Description: Human recombinant full-length AHA1 (GenBank Accession No. NM_012111), with C-terminal His-tag, MW=38 kDa, expressed in a Baculovirus-infected Sf9 cell expression system.

Assay Conditions: The addition of AHA1 increased ATPase activity of HSP90 α from 0.18 to 0.94 pmol/min/ μg . 50 μl reaction mix (100 mM Tris-HCl pH7.4, 20 mM KCl, 6 mM MgCl₂, 200 μM ATP, and 5 μg Hsp90 α) was incubated with Aha1 (0-10 μg) for 3 hrs at 37 $^{\circ}\text{C}$. 100 μl malachite green reagent was added to stop the reaction. Absorbance was measured at 620 nm.

Application: Useful for the study of HSP90 α activity and biochemical function.

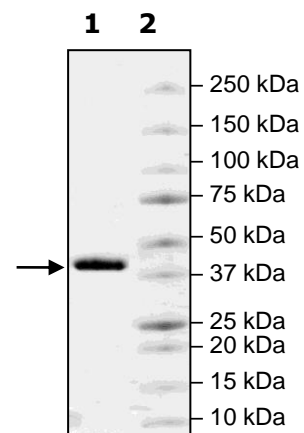
Quality Assurance



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

Lane 1:
2.5 μg Aha1
Lane 2:
Protein Marker

MW: 38 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