

TRAP1 (human), (recombinant) (His- tag)

The Hsp90 family of heat shock proteins represents one of the most abundantly expressed and highly conserved families of cellular chaperones whose expression can be upregulated under conditions of cellular stress. It includes cytoplasmic (Hsp90-alpha/beta), ER (Grp94) and mitochondrial (TRAP1) localized members. Hsp90 is structurally composed of an N-terminal ATP binding domain, a medial substrate-binding region, and a C-terminal dimerization motif. Hsp90 dimers function in cooperation with cochaperones to stabilize a multitude of client proteins substrates. TRAP1 (aka Hsp75) was originally identified as a chaperone binding partner for retinoblastoma protein. TRAP1 was later found by immunofluorescence to be localized to the mitochondria. TRAP1 is significantly more active as an ATPase due in part to the lack of regulatory elements in the C-terminus, but retains the ability to dimerize. However, it does not form stable complexes with typical cochaperones of Hsp90 like p23 and HOP4. TRAP1 has been identified as a protective element for cell survival, and has been suggested as a target for anti-cancer therapeutics.

Citations: 5

[View Online »](#)

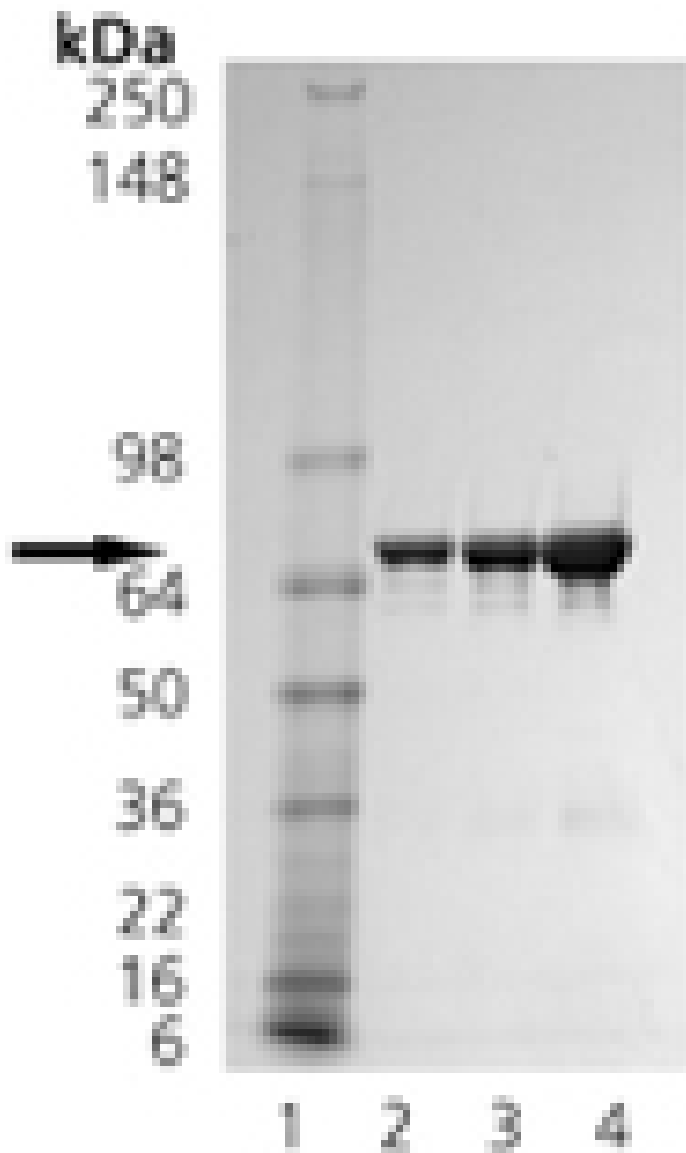
Ordering Information

[Order Online »](#)

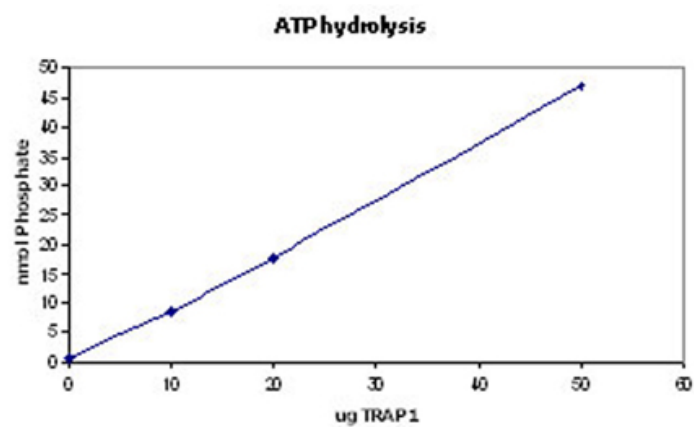
ADI-SPP-848-050	50µg
ADI-SPP-848-200	200µg

Manuals, SDS & CofA

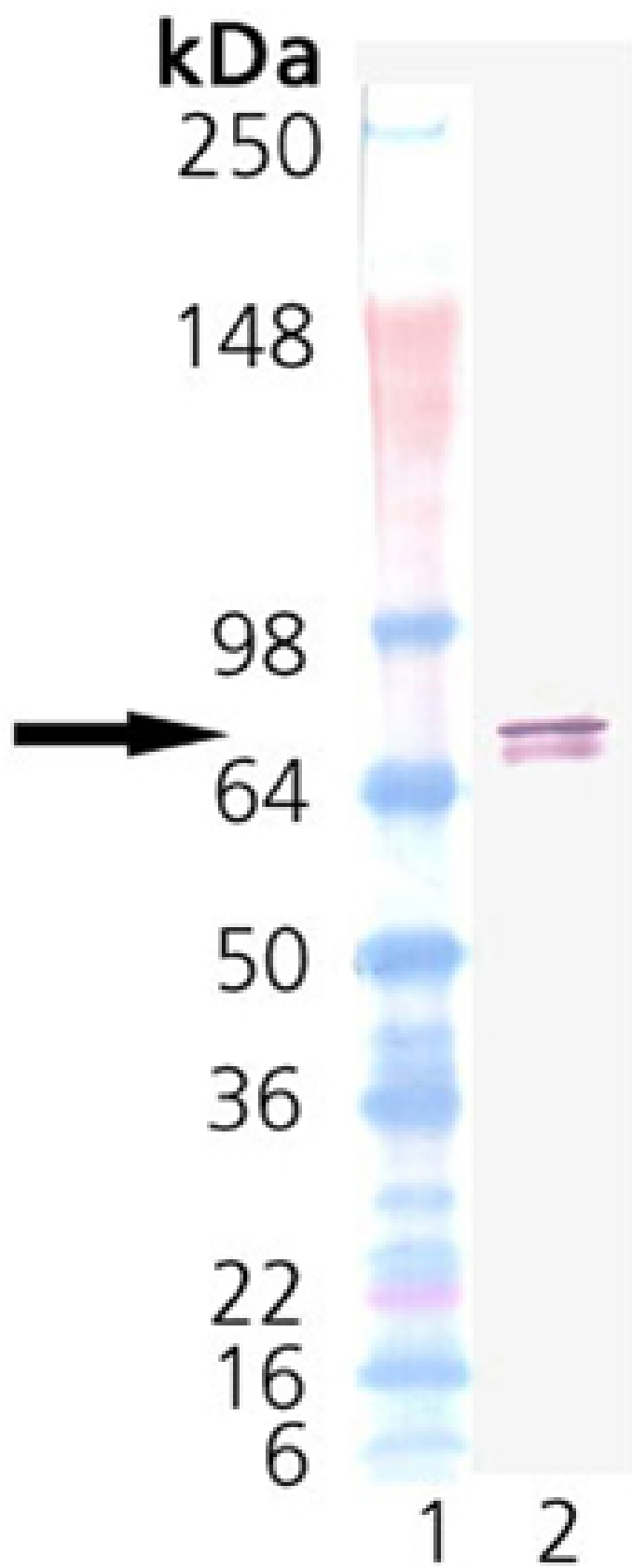
[View Online »](#)



SDS-PAGE analysis: Lane 1: MW marker, Lane 2: 1µg, Lane 3: 2µg, Lane 4: 5µg TRAP1 protein



ATPase Assay: ATP hydrolysis after 1 hour at 37°C detected by malachite green. Quantities determined by comparison to potassium phosphate standard. Approximate activity is 1.2 nmol PO_4 /nmol TRAP1/min.



Western Blot analysis: Lane 1: MWM, Lane 2: TRAP1, probed with TRAP1 mAb.

Handling & Storage

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name HSP75, Mitochondrial HSP90, MthSP90

Application Notes Western blot control.

Formulation Liquid. In 50mM TRIS, pH 7.5, containing 150mM sodium chloride, 1mM DTT, and 1mM EDTA.

MW ~72kDa

Purity ≥95% (SDS-PAGE; Western blot)

Purity Detail Purified by multi-step chromatography.

Source Produced in *E. coli*. Human TRAP1 (aa 60-704) is fused at the N-terminus to a His-tag.

UniProt ID Q12931



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)