

# HSP60 monoclonal antibody (LK-2)

Hsp60 is a member of the chaperonin family of heat shock proteins, with homologs functioning in the cytosol and mitochondria to fold nascent and aggregated proteins. Hsp60 is the eukaryotic homolog of the *E. coli* GroEL protein, and forms a multimeric complex in the mitochondria with Hsp10 (Cpn10) to form a large central cavity in which ATP-dependent protein folding takes place. TRiC/CCT, a eukaryotic relative of Hsp60, is expressed in the cytosol and participates in the folding of actin and tubulin substrates, but lacks any association with an Hsp10-like co-factor.

This antibody is covered by our [Worry-Free Guarantee](#).

Citations: 36

[View Online »](#)

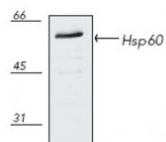
## Ordering Information

[Order Online »](#)

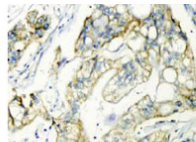
ADI-SPA-807-E	100µg
---------------	-------

## Manuals, SDS & CofA

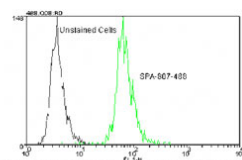
[View Online »](#)



Western blot analysis of pig kidney ESK-4 (heat shocked) cell lysate, probed with Hsp60 mAb (LK-2)



Human colon cancer tissue immunohistochemically stained using Hsp60 Monoclonal Antibody (LK-2)



Flow cytometry analysis of  $10^5$  Jurkat cells using Hsp60 Monoclonal Antibody (LK-2), DyLight™ 488 Conjugate at a concentration of 50  $\mu$ g/ml.

## Handling & Storage

**Long Term Storage**                      -20°C

Shipping Blue Ice

## Regulatory Status RUO - Research Use Only

## Product Details

<b>Alternative Name</b>	Chaperonin 60, CPN60, HspD1, heat shock protein 60
-------------------------	--

<b>Application</b>	ELISA, Flow Cytometry, IHC, IP, WB
--------------------	------------------------------------

<b>Application Notes</b>	Detects a band of ~60kDa by Western blot.
--------------------------	---

**Clone** LK-2

<b>Formulation</b>	Liquid. In PBS, pH 7.2, containing 0.1mM PMSF and 50% glycerol.
--------------------	---

**GenBank ID** M34664

Host Mouse

<b>Immunogen</b>	Recombinant human HSP60.
------------------	--------------------------

**Isotype** IgG1

**Purity Detail** Protein G affinity purified.

<b>Recommendation Dilutions/Conditions</b>	Flow Cytometry (10µg/ml)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
--	---

<b>Source</b>	Purified from ascites.
---------------	------------------------

**Species Reactivity** Bovine, C. elegans, Chicken, Dog, Ehrlichia sennetsu, Guinea pig, Hamster, Human, Monkey, Mouse, Plant, Porcine, Rabbit, Rat

UniProt ID P10809

Worry-free Guarantee

This antibody is covered by our [Worry-Free Guarantee](#)

Last modified: May 29, 2024



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)