# HSP90α polyclonal antibody (PE conjugate)

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, and includes cytoplasmic (Hsp90-α/β), ER (grp94), and mitochondrial (TRAP1) localized members. Structurally, Hsp90 is characterized by an N-terminal ATP-binding domain, a medial substrate-binding domain, and a C-terminal dimerization motif. Hsp90 dimers function in cooperation with cochaperones (e.g. Hsp40, Hsp70, Hop, p23) to stabilize a multitude of client protein substrates, including steroid hormone receptors, protein kinases, and transcription factors. The essential binding and hydrolysis of ATP by Hsp90 is inhibited by ansamycin drugs (e.g. geldanamycin, 17-AAG) which occupy the N-terminal Hsp90 nucleotide-binding pocket. Many Hsp90 client proteins such as erbB2/Her-2, c-raf, bcr-abl, p53, and hTERT, are members of well characterized oncogenic pathways, making Hsp90 inhibitors useful anticancer agents.

This antibody is covered by our Worry-Free Guarantee.

Citations: 2

View Online »

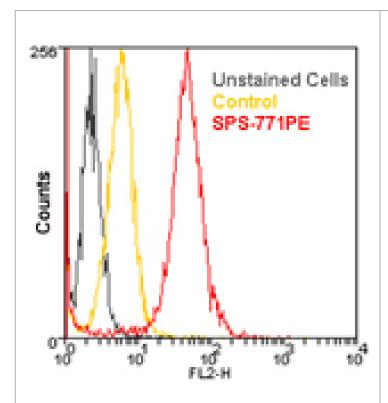
# **Ordering Information**

**Order Online** »

ADI-SPS-771PE-050	50µg
ADI-SPS-771PE-200	200μg

Manuals, SDS & CofA

**View Online** »



Flow cytometry analysis of  $10^6$  Jurkat cells using Hsp90 $\alpha$ , pAb (PE conjugate) at a concentration of 10  $\mu$ g/ml.

## **Handling & Storage**

**Handling** Avoid vigorous centrifugation and vortexing. Do not freeze. Protect from light.

Long Term Storage +4°C

Shipping Blue Ice

# Regulatory Status RUO - Research Use Only

### **Product Details**

Alternative Name HSP86, Heat shock protein 90α

**Application** Flow Cytometry

**Formulation** Liquid. In PBS, pH 7.2, containing 0.09% sodium azide.

GenBank ID X16857

**Host** Rabbit

Immunogen Synthetic peptide corresponding to a portion of mouse

Hsp86.

Purity Detail Protein A affinity purified.

Recommendation Dilutions/Conditions Flow Cytometry (1:100)Suggested dilutions/conditions may

not be available for all applications. Optimal conditions must be determined individually for each application.

Source Purified from rabbit serum.

**Species Reactivity** Beluga, Bovine, Dog, Fish, Guinea pig, Hamster, Human,

Monkey, Mouse, Porcine, Rabbit, Rat, Sheep, Xenopus

UniProt ID P07901

### Worry-free Guarantee This antibody is covered by our Worry-Free Guarantee



European Sales Office ENZO LIFE SCIENCES (ELS) AG Phone: +41 61 926 8989 infoeu@enzolifesciences.com Belgium, The Netherlands & Luxembourg Phone: +32 3 466 0420 infobe@enzolifesciences.com

Phone: +33 472 440 655 infofr@enzolifesciences.com Germany
Phone: +49 7621 5500 526
infode@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
infouk@enzolifesciences.com