# HSP104 polyclonal antibody

This antibody is covered by our Worry-Free Guarantee.

Citations: 2

View Online »

**Ordering Information** 

Order Online »

ALX-210-140-R100

100µl

Manuals, SDS & CofA

**View Online »** 

## **Handling & Storage**

**Handling** Avoid freeze/thaw cycles.

Long Term Storage -20°C

Shipping Blue Ice

# Regulatory Status RUO - Research Use Only

### **Product Details**

Alternative Name Heat shock protein 104

**Application** WB

**Application Notes**Detects a band of ~104kDa (yeast). Does not cross-react

with HSP100 from prokaryotes, nor certain plants or

insects including Drosophila.

Crossreactivity Does not cross-react with HSP100 from prokaryotes, nor

certain plants or insects including Drosophila.

**Formulation** Liquid. Contains 0.05% sodium azide.

**Host** Rabbit

**Immunogen** Synthetic peptide corresponding to aa 894-908

(D894DDNEDSMEIDDDLD908) of C-terminal S. cerevisiae HSP104 (heat shock protein 104).

**Recommendation Dilutions/Conditions**Western Blot (1:5,000)Suggested dilutions/conditions may

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

Species Reactivity Yeast

UniProt ID P31539

### **Worry-free Guarantee**

ENZO LIFE SCIENCES,
European Sales Office
Belgium, The Netherlands
France

INC.
ENZO LIFE SCIENCES
& Luxembourg
Phone: +33 472 440 655

Phone: 800.942.0430
(ELS) AG
Phone: +32 3 466 0420
info

 $\begin{tabular}{lll} \underline{usa@enzolifesciences.com} & & \underline{info} \\ & & \underline{eu@enzolifesciences.com} \end{tabular}$ 

Phone: +41 61 926 8989

infobe@enzolifesciences.com Phone: +49 7621 5500 526 infode@enzolifesciences.com

This antibody is covered by our Worry-Free Guarantee

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

Last modified: May 29, 2024

info-