# CHO host cell proteins polyclonal antibody

Host cell proteins (HCPs) are proteins unique to the host cell used in biologics production and are considered to be a potentially hazardous contaminant when included in the final recombinant product. There is a range from 4,300 genes in *E. coli* to 30,000 genes in CHO or NS0 cell lines that may be expressed along with the desired recombinant product, and at least a small sub-population of these proteins will normally co-purify with the protein of interest. Characterization of these HCPs is important because there are potential safety risks when they are included in a therapeutic product, and the foreign nature of the proteins can elicit an immune response in humans. Antibodies raised against host cell proteins can be used to detect HCPs using Western blot analysis.

This antibody is covered by our Worry-Free Guarantee.

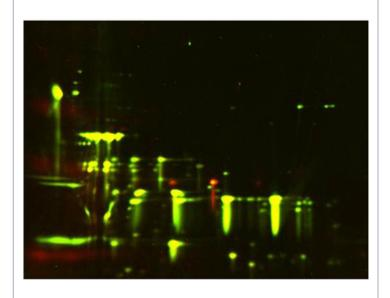
# **Ordering Information**

Order Online »

**ENZ-ABS260-0100** 100μg

Manuals, SDS & CofA

**View Online** »



Western Blot Analysis of CHO HCP, 2-D Gel Separation.

## **Handling & Storage**

**Use/Stability** Stable for one month at 4°C, stable for at least 12 months at -20°C.

**Handling** Avoid freeze/thaw cycles.

Short Term Storage +4°C

Long Term Storage -20°C

Shipping Blue Ice

### Regulatory Status RUO - Research Use Only

### **Product Details**

Alternative Name HCP

**Application** WB

**Formulation** Liquid. In PBS only.

**Host** Rabbit

**Immunogen** CHO host cell proteins.

Purity Detail Protein A affinity purified.

Species Reactivity Hamster

**Specificity** Recognizes CHO host cell proteins.

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 infousa@enzolifesciences.com 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

France
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 info-

uk@enzolifesciences.com