

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](#).

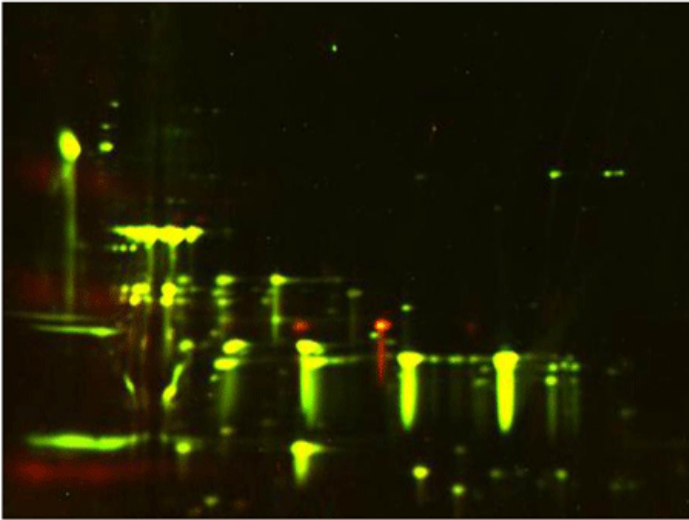
Ordering Information

[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
[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-
fr@enzolifesciences.com](mailto:info-fr@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)