IL-23A polyclonal antibody

This antibody is covered by our Worry-Free Guarantee.

Ordering Information

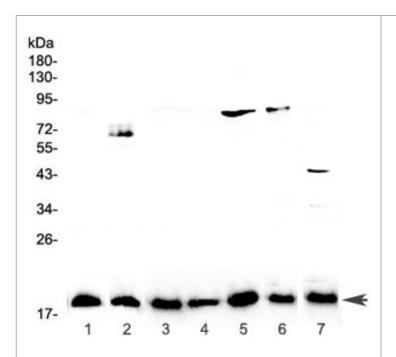
Order Online »

ENZ-ABS898-0100

100µg

Manuals, SDS & CofA

View Online »



Western blot testing of 1) placenta, 2) PANC-1, 3) rat lymph, 4) rat small intestine, 5) rat testis, 6) rat ovary, and 7) mouse testis lysate with IL-23 antibody at 0.5µg/ml. Predicted molecular weight ~21 kDa, observed molecular weight ~19 kDa.

Handling & Storage

Use/Stability After reconstitution, the IL-23 antibody can be stored for up to one month at 4°C. For

long-term, aliquot and store at -20°C.

Handling Avoid freeze/thaw cycles.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Interleukin-23 subunit α, IL-23 p19

Application ELISA, WB

Formulation Lyophilized from 1X PBS containing 2% Trehalose and

0.025% sodium azide.

Host Rabbit

Immunogen Recombinant human IL-23 (aa 20-178).

Isotype IgG

Purity Detail Antigen affinity purified.

Recommendation Dilutions/Conditions ELISA (direct) (0.1-0.5µg/ml)Western blot (0.5-

1μg/ml)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be

determined individually for each application.

Species Reactivity Human, Mouse, Rat

UniProt ID Q9NPF7

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

•

Last modified: July 12, 2024



info-

eu@enzolifesciences.com