

Transferrin polyclonal antibody

Transferrin is a serum glycoprotein involved in iron transport. Transferrin exists in two forms, the iron-free Apotransferrin, which binds two Fe³⁺ ions to generate Ferrotransferrin. Apotransferrin remains tightly bound to its receptor until the complex is recycled to the cell surface, where Apotransferrin is released into the extracellular space to recruit more Fe³⁺ ions.

This antibody is covered by our [Worry-Free Guarantee](#).

Citations: 1

[View Online »](#)

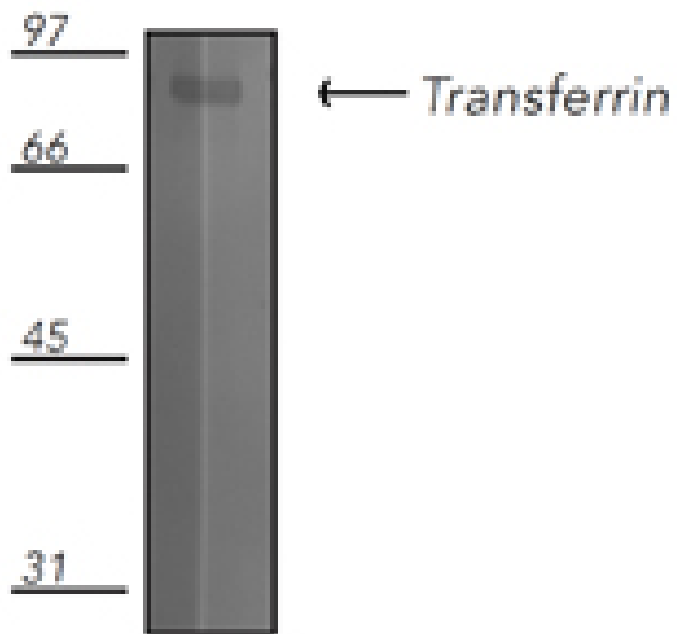
Ordering Information

[Order Online »](#)

ADI-VAP-EN004-D	50µg
ADI-VAP-EN004-F	200µg

Manuals, SDS & CofA

[View Online »](#)



Western blot analysis of rat apotransferrin protein probed with transferrin pAb.

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 TF

Application IP, WB

Application Notes Detects a band of ~76kDa by Western blot.

Formulation Liquid. In PBS containing 50% glycerol and 0.09% sodium azide.

Host Rabbit

Immunogen Human Transferrin.

Purity Detail Protein A affinity purified.

Recommendation Dilutions/Conditions Western Blot (1:2,000, ECL) 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 Dog, Guinea pig, Human, Mouse, Rat

UniProt ID P02787

Worry-free Guarantee This antibody is covered by our [Worry-Free Guarantee](#).



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)