

# Tapasin polyclonal antibody

Tapasin is a 48 kDa glycoprotein which is a member of the immunoglobulin superfamily and is required for efficient peptide binding to TAP (Transporter associated with antigen processing). TAP binds peptides in its cytosolic part and subsequently translocates the peptides into the lumen of the endoplasmic reticulum where assembly of MHC class I and peptide takes place.

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

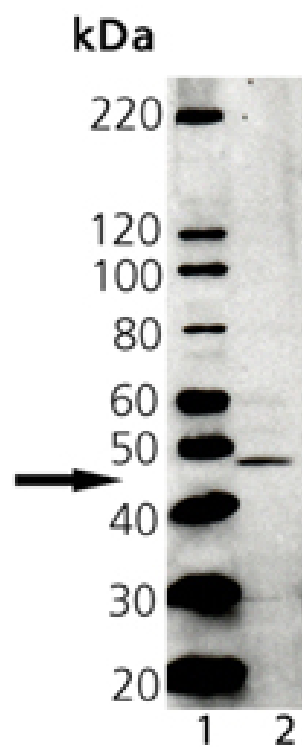
## Ordering Information

[Order Online »](#)

ADI-CSA-630-D	50µg
ADI-CSA-630-F	200µg

## Manuals, SDS & CofA

[View Online »](#)



Western blot analysis of Tapasin, PAb: Lane 1: MW marker, Lane 2: A431 Lysate (Prod No.: ADI-LYC-4100)

# Handling & Storage

Long Term Storage                    -20°C

Shipping                                Blue Ice

## Regulatory Status

RUO - Research Use Only

## Product Details

Alternative Name	TPN, TPSN, TAP binding protein
Application	IP, WB
Application Notes	Detects a band of ~48kDa by Western blot.
Crossreactivity	Weakly cross-reacts with bovine Tapasin.
Formulation	Liquid. In PBS containing 50% glycerol and 0.09% sodium azide.
GenBank ID	AF009510
Host	Rabbit
Immunogen	Synthetic peptide corresponding to aa 2-20 of human tapasin with a cysteine residue added.
Purity Detail	Protein A affinity purified.
Recommendation Dilutions/Conditions	Immunoprecipitation (1:100)Western Blot (1:1,000, colorimetric and 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	Human
UniProt ID	O15533
Worry-free Guarantee	This antibody is covered by our <a href="#">Worry-Free Guarantee</a>



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-  
be@enzolifesciences.com](mailto:info-be@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)