

# Estrogen receptor $\beta$ polyclonal antibody

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

Citations: 1

[View Online »](#)

Manuals, SDS & CofA

[View Online »](#)

## 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	ER $\beta$
Application	IF, IHC (PS), WB
Application Notes	Detects a band of ~55kDa by Western blot.
Crossreactivity	Does not cross-react with ER $\alpha$ .
Formulation	Liquid. In PBS containing 1mg/ml BSA and 0.05% sodium azide.
Host	Rabbit
Immunogen	Synthetic peptide corresponding to aa 467-485 (C467CSTEDSKNKEGSQNLQSQ485) of C-terminal mouse ER $\beta$ (estrogen receptor $\beta$ , isoform $\beta$ -6).
Purity Detail	Epitope-affinity purified.
Recommendation Dilutions/Conditions	Immunofluorescence (1:50-1:200)Immunohistochemistry (paraffin sections, 1-2 $\mu$ g/ml)Western Blot (2 $\mu$ g/ml)
Species Reactivity	Human, Rat
UniProt ID	O08537

## 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-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)

Last modified: May 29, 2024