

# Estrogen receptor $\beta$ polyclonal antibody

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

Citations: 3

[View Online »](#)

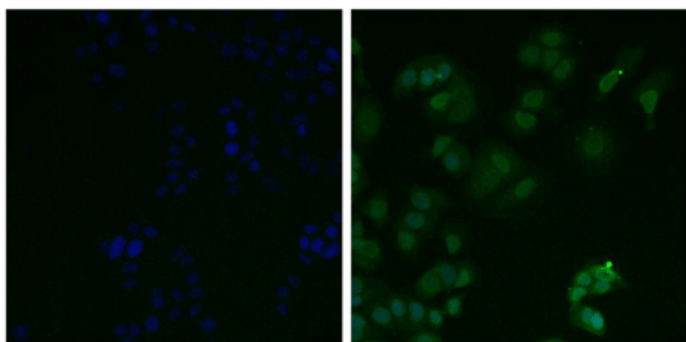
## Ordering Information

[Order Online »](#)

ALX-210-132-C100	100 $\mu$ g
------------------	-------------

Manuals, SDS & CofA

[View Online »](#)



-  
Secondary antibody

ER-beta antibody  
Secondary antibody

Figure 1. Immunofluorescent analysis using ER-beta polyclonal antibody (Prod. No. ALX-210-132) shows nuclear localization of ER-beta in MCF7 cells.

# 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	ERβ
Application	ICC, IF, IHC (FS), IHC (PS), WB
Application Notes	Detects a band of ~55kDa by Western blot.
Crossreactivity	Does not cross-react with ERα.
Formulation	Liquid. In PBS containing 1mg/ml BSA and 0.05% sodium azide.
Host	Rabbit
Immunogen	Synthetic peptide corresponding to aa 512-530 (C512SSTEDSKN KESSQNLQSQ530) of C-terminal rat ERβ (estrogen receptor β).
Purity Detail	Epitope-affinity purified.
Recommendation Dilutions/Conditions	Immunofluorescence (1:100-1:1,000)Immunocytochemistry (1:100-1:1,000)Immunohistochemistry (frozen sections, 1:50)Western Blot (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	Q62986

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)