

# Estrogen receptor monoclonal antibody (h-151)

The Estrogen receptors (ER), mainly ERalpha and ERbeta, are members of the nuclear steroid hormone receptor family and are composed of three domains, an N-terminal domain required for maximal transactivation function (TAF), a DNA binding domain, and a steroid binding domain. Estrogen has pleiotropic effects on the growth, development, and maintenance of diverse tissues.

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

Citations: 5

[View Online »](#)

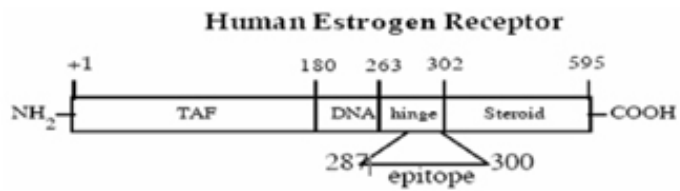
## Ordering Information

[Order Online »](#)

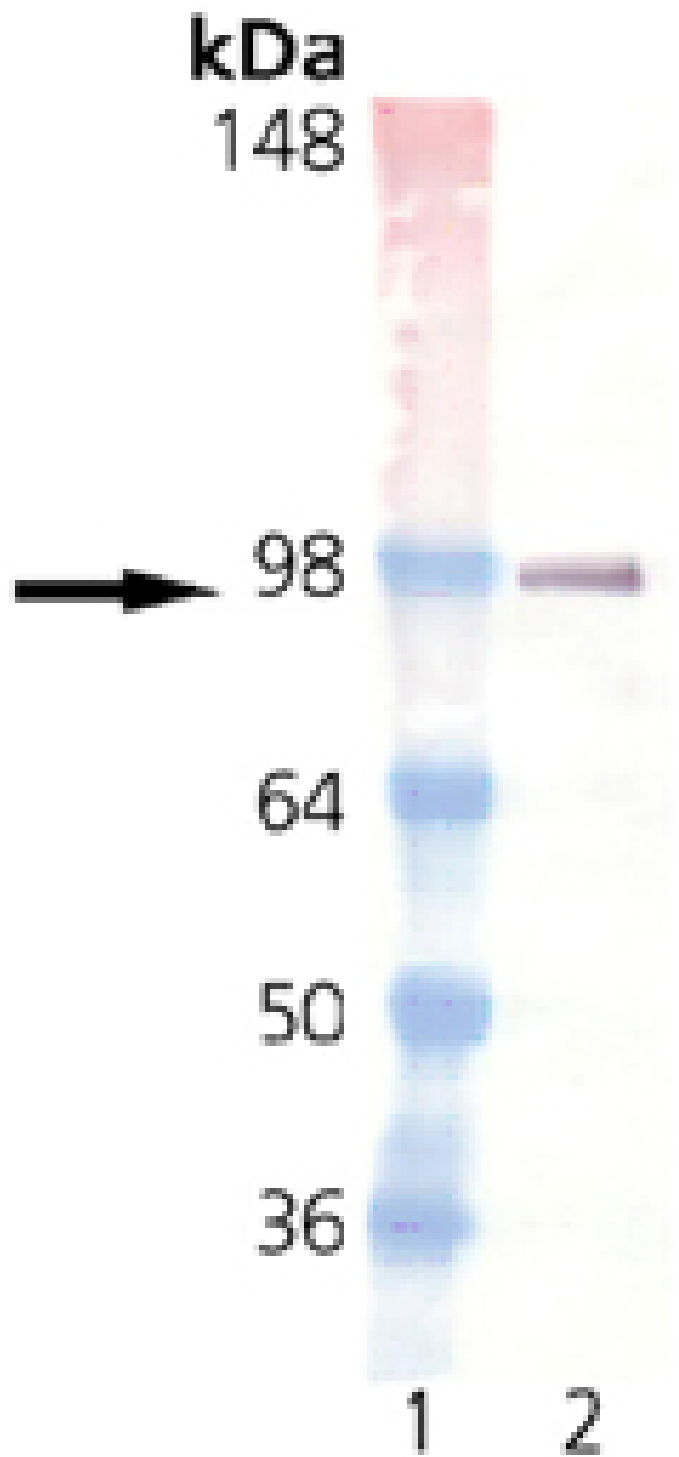
ADI-SRA-1000-D	50µg
ADI-SRA-1000-F	200µg

## Manuals, SDS & CofA

[View Online »](#)



Human Estrogen Receptor domains.



Western blot analysis of ESR1: Lane 1: MWM, Lane 2: Estrogen Receptor 1 (GST-tagged ER1 protein control runs at 92 kDa)

## 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** ESR

**Application** ICC, IP, WB

**Application Notes** Detects a band of ~67kDa by Western blot.

**Clone** h-151

**Formulation** Liquid. In PBS, pH 7.2, containing 50% glycerol and 0.09% sodium azide.

**GenBank ID** M12674

**Host** Mouse

**Immunogen** Synthetic peptide corresponding to the sequence of human estrogen receptor, conjugated to KLH.

**Isotype** IgG1

**Purity Detail** Protein G affinity purified.

**Recommendation Dilutions/Conditions** Immunoprecipitation (15µg/ml)Western Blot (1:250)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.

**Source** Purified from mouse ascites.

**Species Reactivity** Human, Rat

**UniProt ID** P03372

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)