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 »

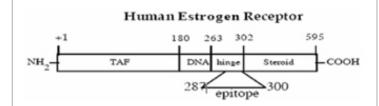
Ordering Information

Order Online »

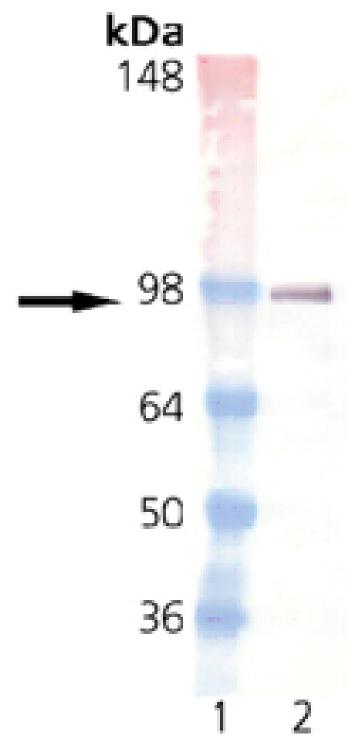
ADI-SRA-1000-D	50µg
ADI-SRA-1000-F	200μg

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 NotesDetects 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.

lsotype lgG1

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

P03372

Worry-free Guarantee

This antibody is covered by our Worry-Free Guarantee

Last modified: May 29, 2024



info-

eu@enzolifesciences.com