

Estrogen receptor α monoclonal antibody (C-542)

The Estrogen receptors (ER), mainly ER α and ER β , 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: 8

[View Online »](#)

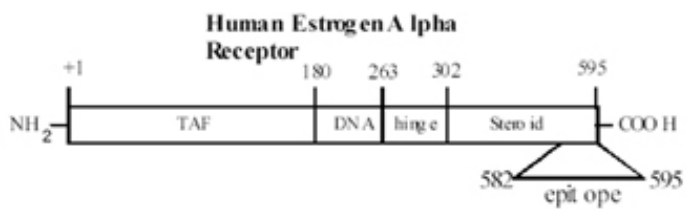
Ordering Information

[Order Online »](#)

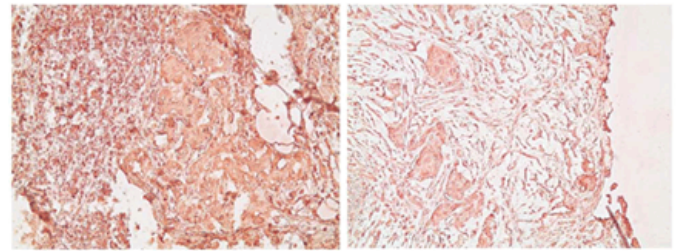
ADI-SRA-1010-D	50 μ g
ADI-SRA-1010-F	200 μ g

Manuals, SDS & CofA

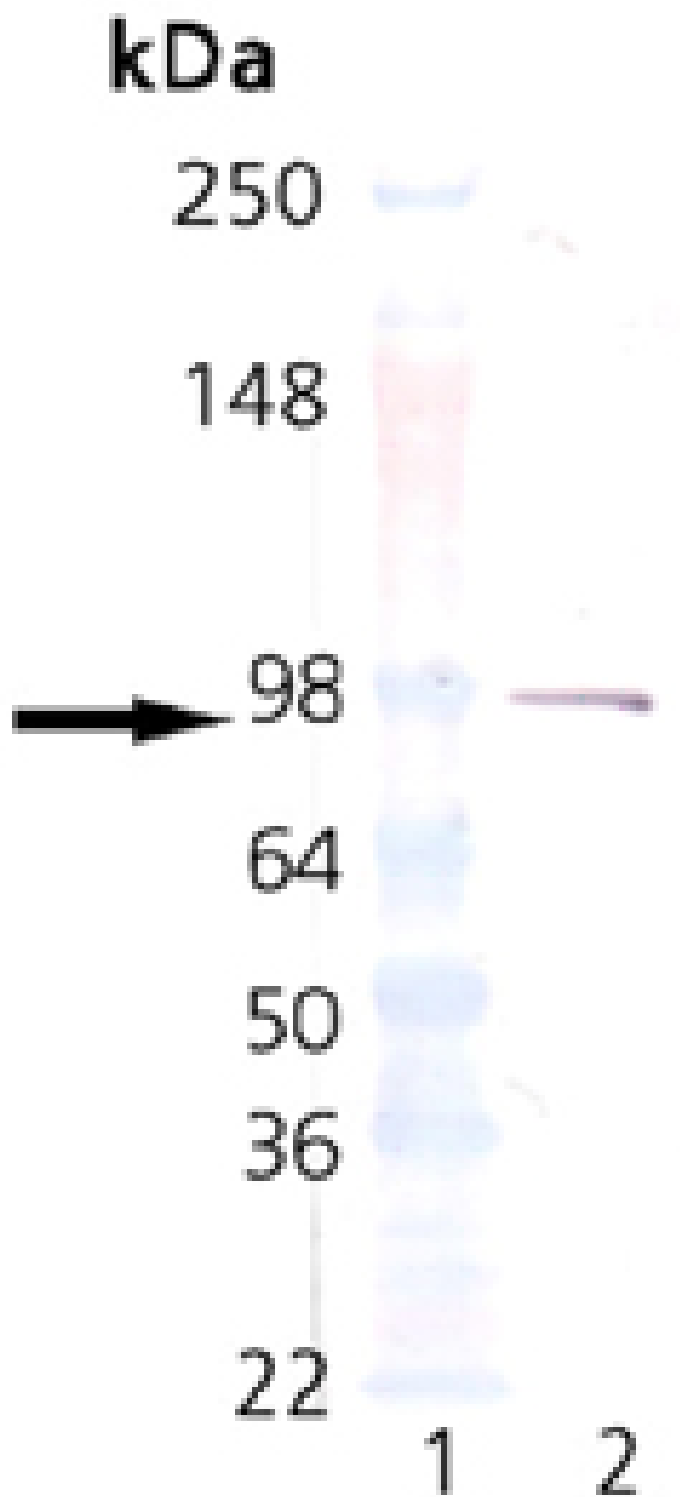
[View Online »](#)



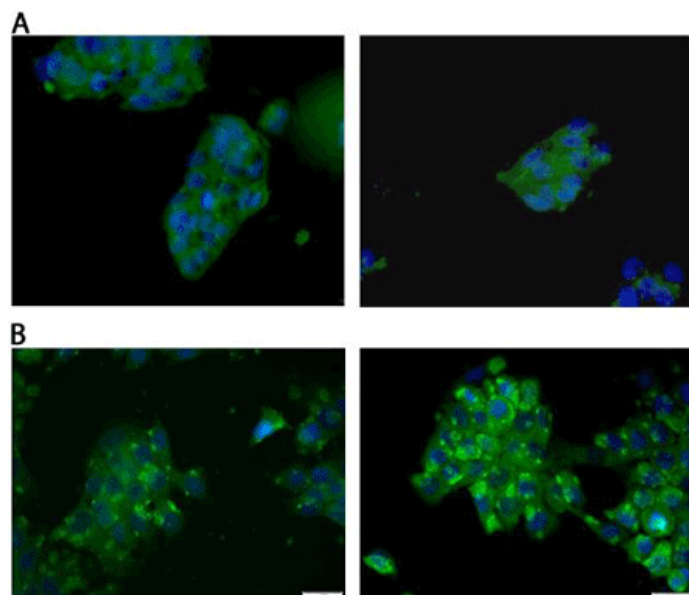
Human Estrogen Alpha Receptor domains.



Immunohistochemistry analysis of human Breast tumor tissue stained with Estrogen receptor α monoclonal antibody (C-542), (Prod. No. ADI-SRA-1010), at 20 μ g/ml.



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



Immunocytochemistry staining of HepG2 cells (Panel A) and MCF-7 (Panel B) with Estrogen receptor α monoclonal antibody (C-542) (Prod. No. ADI-SRA-1010) at 10 μ g/ml.

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, IHC, IP, WB
Application Notes	Detects a band of 66.3kDa by Western blot.
Clone	C-542
Formulation	Liquid. In PBS containing 50% glycerol and 0.09% sodium azide.
GenBank ID	M12674
Host	Mouse
Immunogen	Synthetic peptide corresponding to the sequence near the C-terminus of human Estrogen Receptor α .
Isotype	IgG1
Purity Detail	Protein G affinity purified.
Recommendation Dilutions/Conditions	Western Blot (1:500, colorimetric)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Source	Purified from ascites.
Species Reactivity	Hamster, Human, Mouse, Rat

UniProt ID

P03372

Worry-free Guarantee

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

Last modified: May 29, 2024



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)