

Sox2 recombinant monoclonal antibody (017)

The anti-Sox2 antibody is designed for the detection of human Sox2 by immunohistochemistry.

Sox2 (SRY-box transcription factor 2) is a member of the SOX family of transcription factors characterized by the presence of a high-mobility group (HMG) DNA-binding domain. Sox2 is critically important for the maintenance of pluripotency in embryonic stem cells by regulating genes involved in self-renewal and differentiation. It forms complexes with other transcription factors like Oct4 and Nanog to activate or repress target gene expression.

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

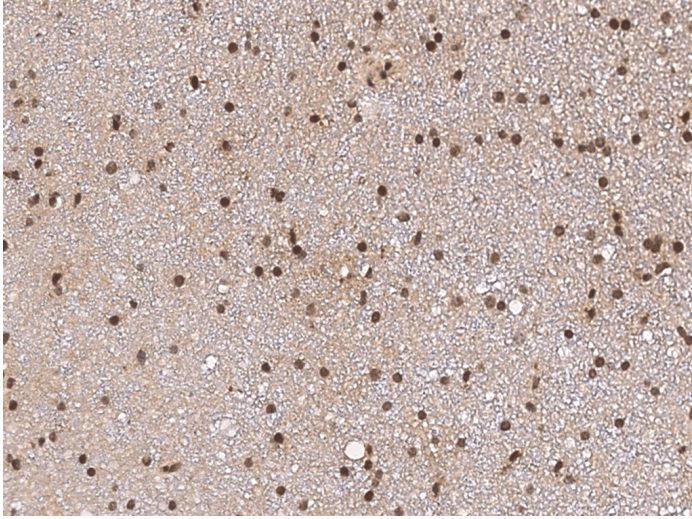
Ordering Information

[Order Online »](#)

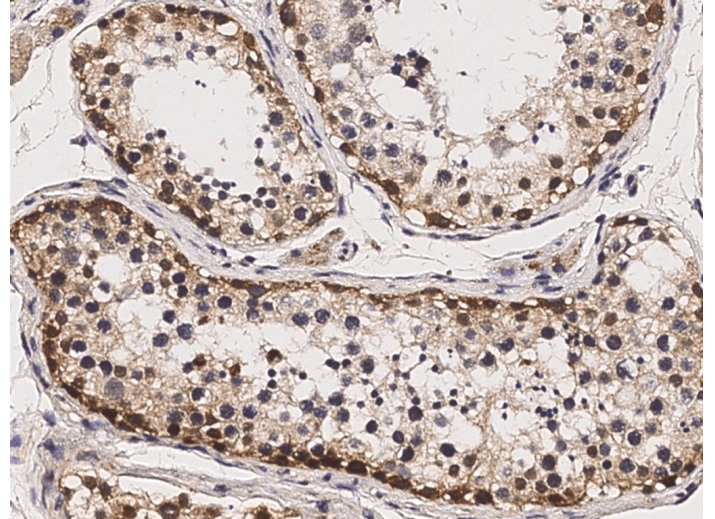
ENZ-ABS1042-0100	100µl
------------------	-------

Manuals, SDS & CofA

[View Online »](#)



Immunohistochemical staining of human SOX2 in human brain with rabbit monoclonal antibody (1:300, formalin-fixed paraffin embedded sections).



Immunohistochemical staining of human SOX2 in human testis with rabbit monoclonal antibody (1:300, formalin-fixed paraffin embedded sections).

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 SRY-box transcription factor 2

Application IHC (PS)

Clone 017

Formulation Liquid. 0.2 µm filtered solution in PBS. Carrier-free. Preservative-free.

Host Rabbit

Immunogen Synthetic peptide corresponding to the N-terminus of the human Sox2.

Isotype IgG

Purity Detail Protein A purified.

Recommendation Immunohistochemistry (paraffin sections) (1:100-1:500)

Dilutions/Conditions
Suggested dilutions/conditions may not be available for all applications.
Optimal conditions must be determined individually for each application.

Species Reactivity Human

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-
fr@enzolifesciences.com](mailto:info-fr@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)