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

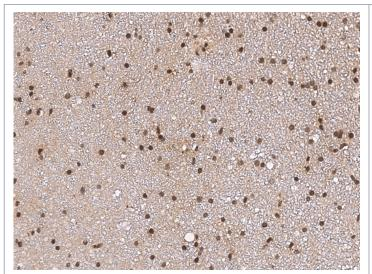
Order Online »

ENZ-ABS1042-0100 10

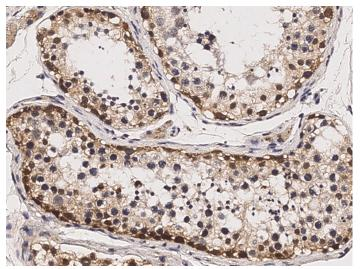
100µl

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)

017 Clone

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

Preservative-free.

Rabbit Host

Synthetic peptide corresponding to the N-terminus of the **Immunogen**

human Sox2.

IgG Isotype

Purity Detail Protein A purified.

Recommendation Dilutions/Conditions Immunohistochemistry (paraffin sections) (1:100-1:500)

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, Phone: 800.942.0430 info-

ENZO LIFE SCIENCES (ELS) AG Phone: +41 61 926 8989 usa@enzolifesciences.com

eu@enzolifesciences.com

European Sales Office

Belgium, The Netherlands & Luxembourg Phone: +32 3 466 0420 infobe@enzolifesciences.com

France Phone: +33 472 440 655 fr@enzolifesciences.com

Phone: +49 7621 5500 526 de@enzolifesciences.com

UK & Ireland Phone (UK customers): 0845 601 1488 Phone: +44 1392 825900 uk@enzolifesciences.com