

HEXB polyclonal antibody

The anti-HEXB antibody is designed to detect human Hexosaminidase subunit beta (HEXB), a critical lysosomal enzyme subunit involved in the degradation of GM2 gangliosides and other glycosaminoglycans. This antibody is optimized for use in immunohistochemistry (IHC) and Western blotting (WB) applications.

HEXB is one of the subunits of the enzyme beta-hexosaminidase, a lysosomal hydrolase responsible for degrading GM2 gangliosides and other glycosaminoglycans. Along with the alpha subunit (encoded by HEXA), HEXB forms functional isoenzymes (HexA and HexB) that play essential roles in lysosomal catabolism. HEXB is expressed in various tissues, with particularly high activity in the central nervous system. Mutations in the HEXB gene are associated with Sandhoff disease, a lysosomal storage disorder characterized by the accumulation of GM2 gangliosides, leading to progressive neurodegeneration. The study of HEXB is critical for understanding the molecular mechanisms of lysosomal function, glycosaminoglycan metabolism, and their dysregulation in diseases such as Sandhoff disease and other neurodegenerative conditions.

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

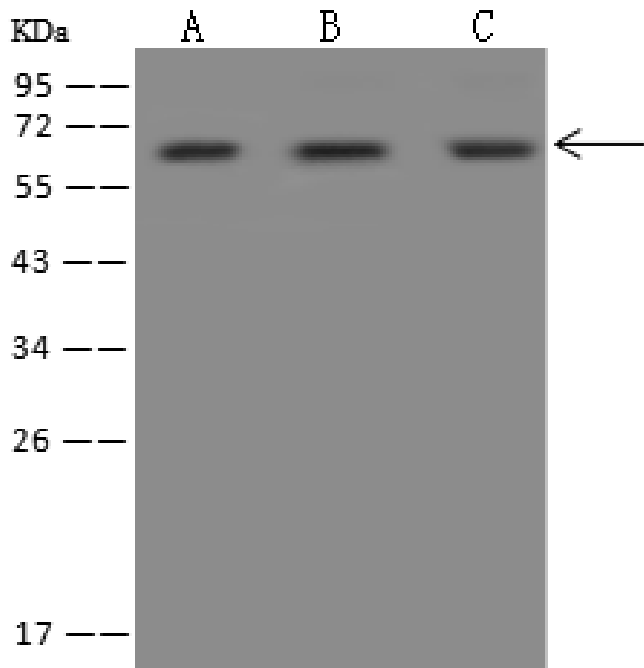
Ordering Information

[Order Online »](#)

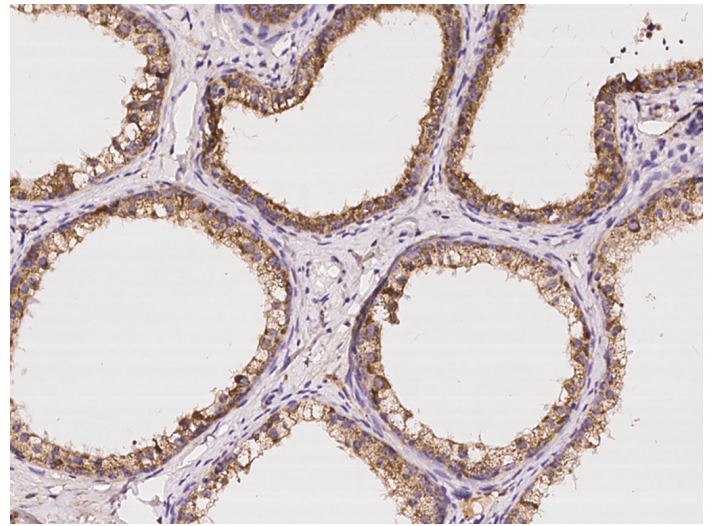
ENZ-ABS1054-0100	100µl
------------------	-------

Manuals, SDS & CofA

[View Online »](#)



Anti-HEXB rabbit polyclonal antibody at 1:500 dilution
Lane A: Jurkat Whole Cell Lysate. Lane B: HeLa Whole Cell Lysate. Lane C: HepG2 Whole Cell Lysate
Lysates/proteins at 30 µg per lane. Secondary ab: Goat Anti-Rabbit IgG (H+L)/HRP at 1/10000 dilution.
Predicted band size: 63 kDa



Immunohistochemical staining of human HEXB in human epididymis with rabbit polyclonal antibody at 1:100 dilution, 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 Beta-hexosaminidase subunit beta

Application IHC (PS), WB

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

Host Rabbit

Immunogen Recombinant human HEXB fragment.

Isotype IgG

Purity Detail Protein A and antigen affinity purified.

Recommendation Dilutions/Conditions Western Blot (1:500-1:2'000)
Immunohistochemistry (paraffin sections) (1:50-1:200)

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,
Last modified January 9, 2025
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

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

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

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

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