

# Hexokinase 3 monoclonal antibody (AT4C12)

Hexokinase is the first enzyme in the glycolytic pathway, catalyzing the transfer of a phosphoryl group from ATP to glucose to form glucose-6-phosphate and ADP. In mammals, four distinct enzymes - types 1 to 4 hexokinases - have been identified. The enzyme is found in most cells, but there is tissue specificity for the particular type of hexokinase. Hexokinase 3 lacks the hydrophobic N-terminal sequence critical for targeting to mitochondria. Hexokinase 3 may have anabolic functions, providing H6P for glycogen or lipid synthesis.

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

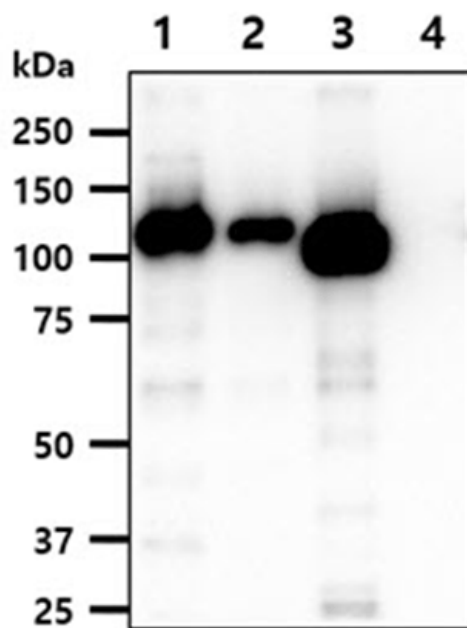
## Ordering Information

[Order Online »](#)

ENZ-ABS757-0100	100µl
-----------------	-------

## Manuals, SDS & CofA

[View Online »](#)



The recombinant protein (20ng) of Hexokinase four isoforms were resolved by SDS-PAGE, transferred to PVDF membrane and probed with Hexokinase 3 antibody (1:1000). Proteins were visualized using a goat anti-mouse secondary antibody conjugated to HRP and an ECL detection system. Lane 1: Recombinant human Hexokinase 1, Lane 2: Recombinant human Hexokinase 2, Lane 3: Recombinant human Hexokinase 3, Lane 4: Recombinant human Hexokinase 4.

## 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** HK3, Hexokinase type III, Hexokinase-C

**Application** ELISA, WB

**Clone** AT4C12

**Formulation** Liquid. In PBS, pH 7.4, containing 0.02% sodium azide and 10% glycerol.

**Host** Mouse

**Immunogen** Recombinant human Hexokinase 3 (aa 1-923) purified from E. coli.

**Isotype** IgG1κ

**Purity Detail** Protein A affinity purified

**Recommendation Dilutions/Conditions** Western Blot (1:2,000). Suggested dilutions/conditions may not be available for all applications. Optimal conditions must be determined individually for each application.

**Species Reactivity** Human

**UniProt ID** P52790

**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)