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. Hexpkinase 3 may have anabolic functions, providing H6P for glycogen or lipid synthesis.

This antibody is covered by our Worry-Free Guarantee.

Ordering Information

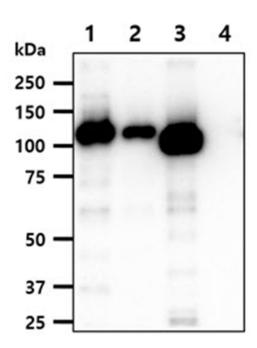
Order Online »

ENZ-ABS757-0100

100µl

Manuals, SDS & CofA

View Online »



The recombianat protein (20ng) of Hexokinase four isoform 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 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.

lgG1k

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



European Sales Office ENZO LIFE SCIENCES (ELS) AG Phone: +41 61 926 8989 infoeu@enzolifesciences.com Belgium, The Netherlands & Luxembourg Phone: +32 3 466 0420 infobe@enzolifesciences.com

Phone: +33 472 440 655 infofr@enzolifesciences.com Phone: +49 7621 5500 526 infode@enzolifesciences.com UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
infouk@enzolifesciences.com