

BEX5 Polyclonal Antibody

(Catalog # A65114)

Background

extracellular exosome

Description

BEX5 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

Liquid. 0.03% Proclin 300, 50% Glycerol, 0.01M PBS, PH 7.4.

Specificity

Human

Isotype

IgG

Uniprot ID

Q5H9J7

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Protein BEX5 protein (1-111AA)

Storage

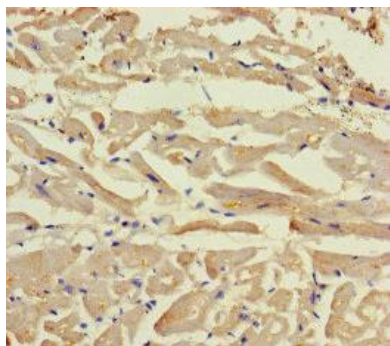
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

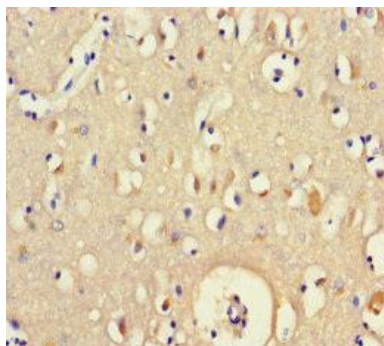
Protein BEX5, Brain-expressed X-linked protein 5, NGFRAP1-like protein 1, Nerve growth factor receptor-associated protein 2, BEX5, NADE2, NGFRAP1L1

Application

ELISA, IHC; Recommended dilution: IHC:1:20-1:200



Immunohistochemistry of paraffin-embedded human heart tissue using BEX5 Polyclonal Antibody at dilution of 1:100



Immunohistochemistry of paraffin-embedded human brain tissue using BEX5 Polyclonal Antibody at dilution of 1:100