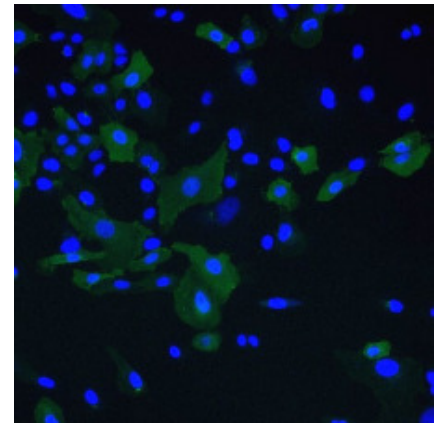


Mouse monoclonal antibody to human Artemin

326-100

Name	Mouse monoclonal antibody to human Artemin
Immunogen	Human Artemin
Immunogen Description	Recombinant human Artemin protein produced using CHO-based Icosagen Cell factory Ltd. proprietary suspension cell line. Immunogen is purified from cell culture supernatant
Uniprot ID	Q5T4W7
Alternative Names	Enovin, Neublastin
Clonality	Mouse monoclonal
Clone	2G7
Class	IgG1
Reactivity	Human Artemin
Application	ELISA, IF, IHC
Protocol	ELISA 0.2 to 1 µg/ml; IF 1 to 20 µg/ml; IHC 5 µg/ml. Conformational antibody, not suitable for Western Blot application. Monoclonal antibody working titer has to be established practically for each particular antigen and assay format
Purification	Protein G purification
Buffer	PBS pH 7.4, with 0.1% sodium azide
Unit Size	100 µg
Related Products	Monoclonal and polyclonal antibodies

100µg



B

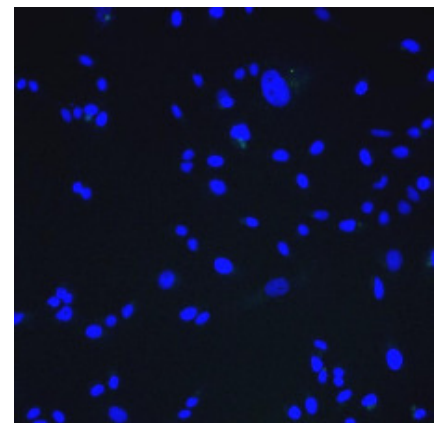


Figure 1. Immunofluorescence analysis of GFRA3-Ret51 receptor complex bound Artemin. Human U2OS cells were transfected with GFRa3-Ret51 expression vector followed by treatment with 100 nM Artemin. Artemin bound to the GFRa3-Ret51 receptor complex was detected by anti-Artemin antibody clone 2G7. Goat anti-mouse Dylight 488 conjugated antibody (dilution 1:1000) was used as secondary antibody. For nuclear staining DAPI was used. ArrayScan VTI platform (Thermo Scientific) was used

and ELISA kit to human Artemin. For more information visit www.icosagen.com/antibodies or <http://icosagen.com/products/elisa-kits>

Shipping

This product is shipped in non-frozen liquid form in ambient conditions

Storage

Store at $-20...-70^{\circ}\text{C}$ upon receipt. Divide antibody into aliquots prior usage. Avoid multiple freeze-thaw cycles

Background

Artemin supports the survival of sensory and sympathetic peripheral neurons in culture and also supports the survival of dopaminergic neurons of the ventral mid-brain. Strong attractant of gut hematopoietic cells thus promoting the formation Peyer's patch-like structures, a major component of the gut-associated lymphoid tissue. Ligand for the GFR-alpha-3-RET receptor complex but can also activate the GFR-alpha-1-RET receptor complex (UniProt)

Limitations

This product is for research use only

TECHNICAL ASSISTANCE

Please refer any technical questions to technical.support@icosagen.com

For Research Use Only

distributed in the US/Canada by:
EAGLE BIOSCIENCES, INC.
20A NW Blvd, Suite 112 Nashua, NH 03063
Phone: 617-419-2019 FAX: 617-419-1110
www.EagleBio.com • info@eaglebio.com



for image acquisition (10x objective). Composite picture was generated using pseudocolors green for Artemin and blue for nuclei. A. Artemin bound to the GFR α 3-Ret51 receptor complex; B. Negative control - GFR α 3-Ret51 receptor expressing U2OS cells

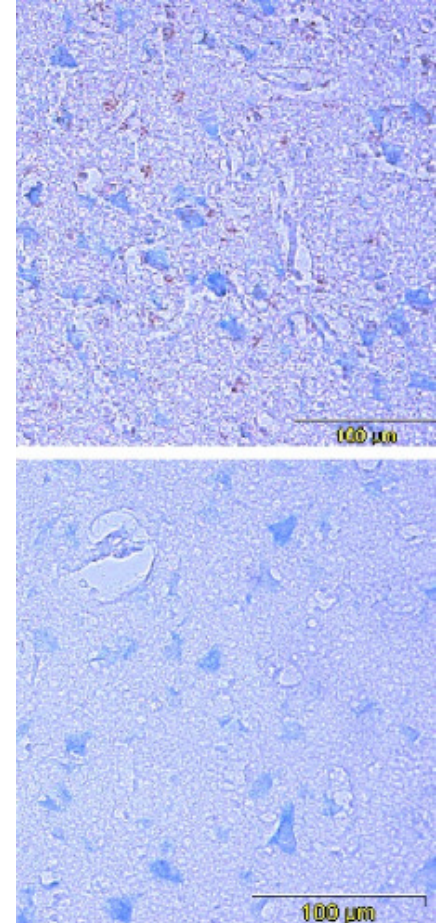


Figure 2.
Immunohistochemistry testing of anti-ARTN monoclonal antibody 2G7.
Analysis was performed using FFPE human cerebral cortex tissue sections from Alzheimer's disease patients. Tissue sections were boiled with sodium citrate buffer (pH 6) for antigen retrieval. Incubation with primary antibody at 5 $\mu\text{g}/\text{ml}$ was performed overnight at 4 $^{\circ}\text{C}$. DAB was used for visualization. Sections were counterstained with toluidine blue.