



DATA SHEET

Mouse monoclonal antibody to human Artemin

Catalog no: 325-100**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:** 1F3**Class:** IgG1**Reactivity:** Human Artemin**Application:** ELISA, IF**Protocol:** ELISA 1:1000 to 1:10 000; IF 1:50 – 1:3000. **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:** IgG fraction purified with Protein G affinity chromatography**Buffer:** PBS pH 7.4, with 0.1% sodium azide.**Concentration:** 1 mg/ml**Unit size:** 100 µg**Related products:** Monoclonal and polyclonal antibodies and ELISA kit to human Artemin.
For more information visit<http://icosagen.com/products/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°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)**References:** -



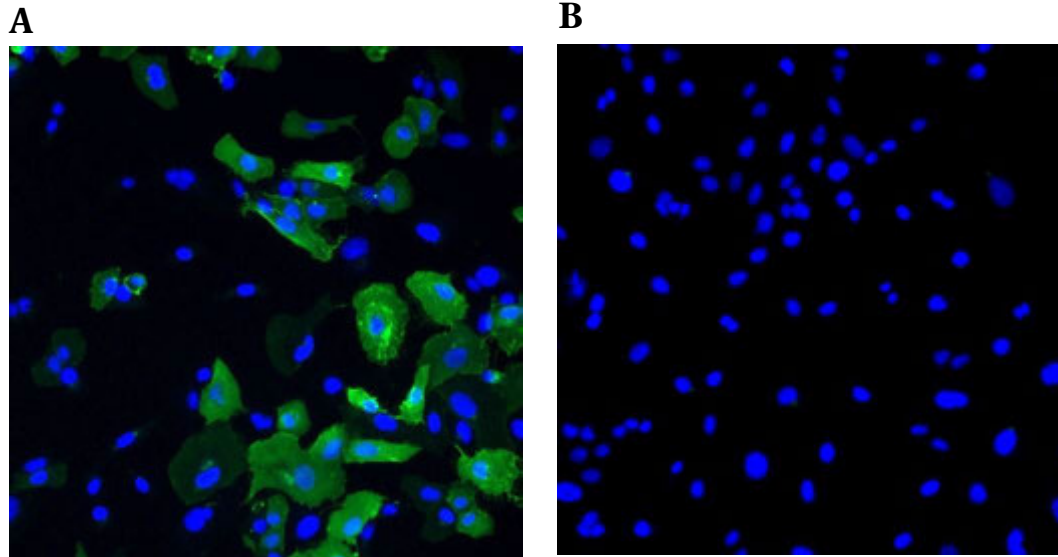


Figure 1. Immunofluorescence analysis of GFRA3-Ret51 receptor complex bound Artemin. Human U2OS cells were transfected with GFRA3-Ret51 expression vector followed by treatment with 100 nM Artemin. Artemin bound to the GFRA3-Ret51 receptor complex was detected by anti-Artemin antibody clone 1F3. Goat 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 for image acquisition (10x objective). Composite picture was generated using pseudocolors green for Artemin and blue for nuclei. A. Artemin bound to the GFRA3-Ret51 receptor complex; B. Negative control - GFRA3-Ret51 receptor expressing U2OS cells.

This product is 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



EAGLE
BIOSCIENCES



Icosagen AS

Eerika tee 1, Õssu küla, Ülenurme vald, Tartumaa 61713, Estonia

Phone +372 737 7070, Fax +372 737 7077, info@icosagen.com

www.icosagen.com