



## DATA SHEET

## Mouse monoclonal antibody to human GDNF

**Catalog no:** 315-100

**Immunogen:** Human Glial cell line-derived neurotrophic factor (GDNF)

**Immunogen description:** Recombinant human GDNF protein produced using *E. coli* expression system

**Uniprot ID:** P39905

**Alternative names:** Astrocyte-derived trophic factor (ATF)

**Clonality:** Mouse monoclonal

**Clone:** 3C1

**Class:** murine IgG1

**Reactivity:** Human GDNF

**Application:** ELISA, WB, IF, IHC

**Protocol:** ELISA 50 to 100 ng/ml; Western immunoblotting 2 to 5 µg/ml ; IF 2 to 10 µg/ml; IHC 5 µg/ml.

Monoclonal antibody working amount 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.

**Unit size:** 100 µg

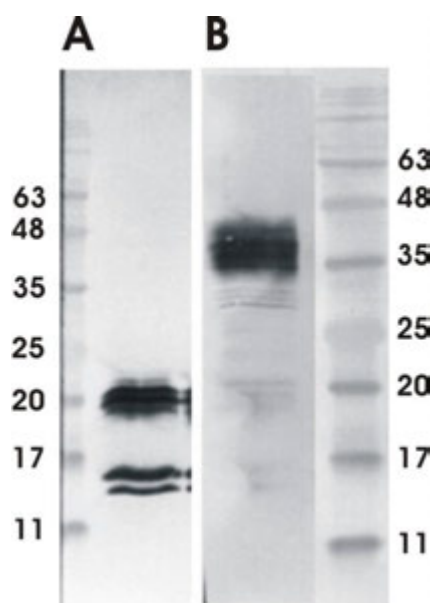
**Related products:** Monoclonal and polyclonal antibodies to human GDNF. For more information visit [www.icosagen.com/antibodies](http://www.icosagen.com/antibodies)

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

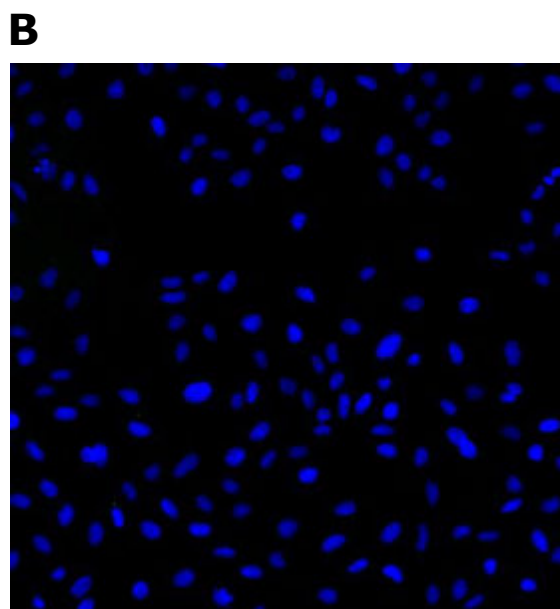
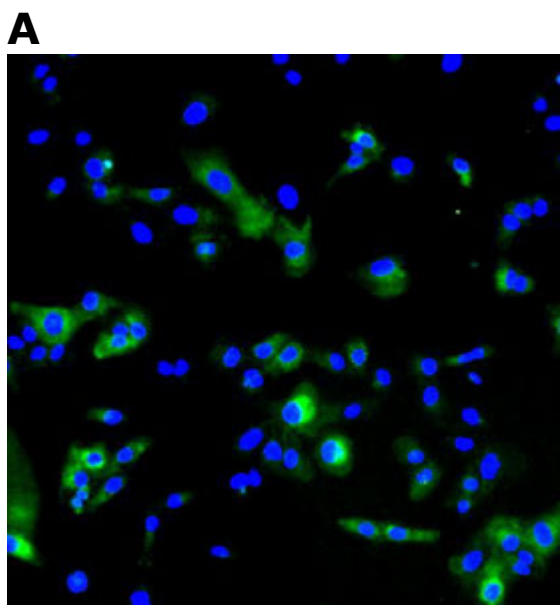
**Storage:** Store at -20...-70°C upon receipt. Divide antibody into aliquots prior usage. Avoid multiple freeze-thaw cycles.

**Background:** GDNF is a neurotrophic factor that enhances survival and morphological differentiation of dopaminergic neurons and increases their high-affinity dopamine uptake. Ligand for the GFR-alpha-3-RET receptor complex but can also activate the GFR-alpha-1-RET receptor complex.

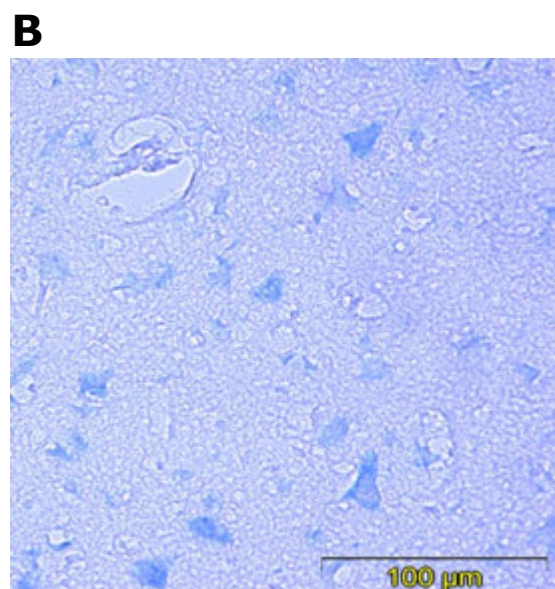
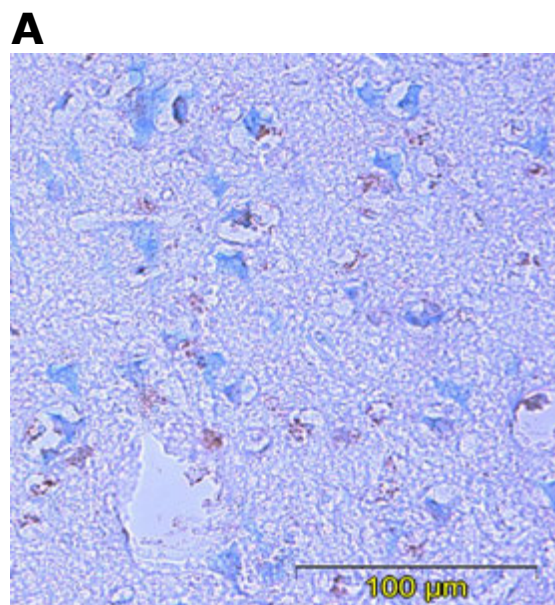
**References:** -



**Figure 1. Western Blot testing of GDNF monoclonal antibody 3C1.** Analysis was performed with antiGDNF monoclonal antibody 3C1 HRP conjugate. 15 µl of CHOEBNALT85 GDNF producing cell culture supernatant was loaded per lane. Analysis was performed in reducing (A) and non-reducing (B) conditions.



**Figure 2. Immunofluorescence detection of human GDNF expressed in U2OS cells by anti-GDNF monoclonal antibody 3C1.** Anti-GDNF antibody 3C1 concentration in IF experiment was 10 µg/ml. Goat anti-mouse AlexaFluor488 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 GDNF specific signal and blue for nuclei. A. GDNF-expressing U2OS cells; B. Negative control (non-transfected U2OS cells)



**Figure 3. Immunohistochemistry testing of anti-GDNF monoclonal antibody 3C1.** Analysis was performed using formalin-fixed paraffin-embedded 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 µg/ml was performed overnight at 4°C. DAKO EnVision™ Detection System, Peroxidase/DAB was used for visualization. Sections were counterstained with toluidine blue and mounted with Eukitt mounting medium. A. GDNF staining by monoclonal antibody 3C1; B. Negative staining without primary antibody

**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](http://www.EagleBio.com) • [info@eaglebio.com](mailto:info@eaglebio.com)

