



DATA SHEET

Mouse monoclonal antibody to human CDNF

Catalog no: 301-100

Immunogen: human CDNF

Immunogen description: Recombinant human CDNF protein produced using CHO-based Icosagen Cell factory Ltd. proprietary suspension cell line. Purified from cell culture supernatant.

Uniprot ID: Q49AH0

Alternative names: ARMETL1

Clonality: Mouse monoclonal

Clone: 7D6

Class: IgG1

Reactivity: Human, mouse; others not tested

Application: ELISA, WB, IF

Protocol:

ELISA - 1:5000 to 1:10 000;

WB - 1:1000 to 1:4000;

IF - 1:100 to 1:3000;

Monoclonal antibody working dilution has to be established practically for each particular antigen and assay format

Purification: purified by protein G affinity chromatography

Buffer: PBS pH 7.4, with 0.1% sodium azide

Concentration: 1 mg/ml

Unit size: 100 µg

Related products: Mono- and polyclonal antibodies to human CDNF. For more details 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 as product degradation may result.

Background: CDNF is a trophic factor for midbrain dopamine neurons *in vivo*. It prevents the 6-OHDA- (Lindholm et al. 20007; Voutilainen et al., 2011) and MPTP-induced degeneration (Airavaara et al., 2012) of dopamine neurons in rodent models of Parkinson's disease. When administered after 6-OHDA or MPTP -lesioning it restores the dopaminergic function and prevents degeneration of dopamine neurons in substantia nigra pars compacta.

References: -

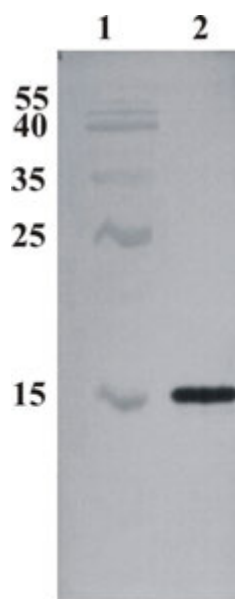


Figure 1. Western Blot testing of anti-CDNF monoclonal antibody (7D6) at 0.25 µg/ml. **Line 1.** PageRuler Prestained Protein Ladder (#SM0671 Fermentas); **Line 2.** Recombinant human CDNF expressed into the supernatant of CHO cell culture medium.



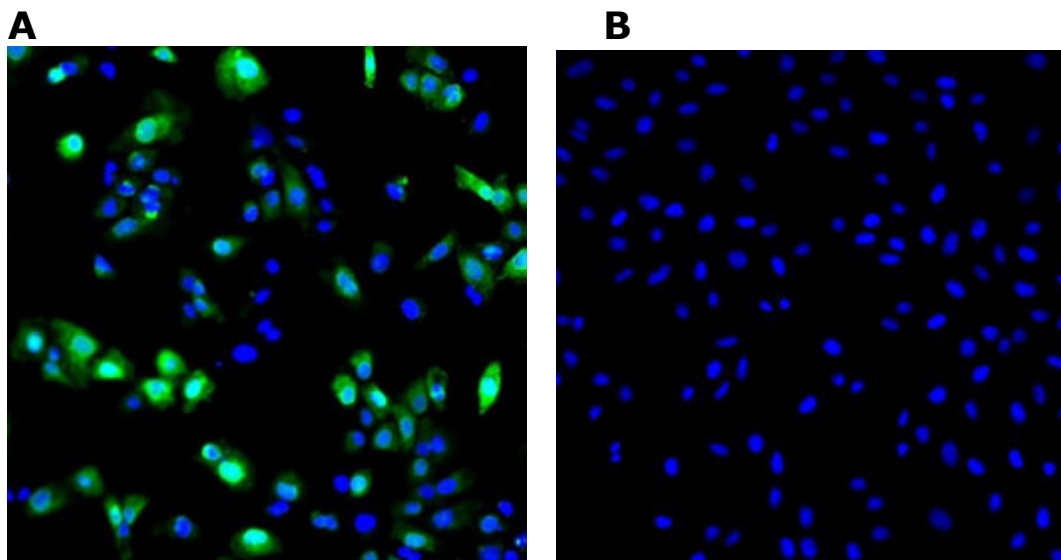


Figure 2. Immunofluorescence detection of human CDNF expressed in U2OS cells. CDNF was visualized using anti-CDNF antibody clone 7D6 at 1 µg/ml. Goat ant-mouse AlexaFluor488 was used as secondary antibody. For nuclear staining DAPI was used. ArrayScan VTI platform (Thermo Scientific) was used for image acquisition (10x objective). Composite picture was generated using pseudocolors green for CDNF specific signal and blue for nuclei. **A.** CDNF-expressing U2OS cells; **B.** Negative control.

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

