

Huntingtin monoclonal antibody - Classic

Other names: HTT, HD, IT15

Cat. No. C15200226

Type: Monoclonal

Isotype: IgG1k

Source: Mouse

Lot #: 001

Size: 50 µg/ 50 µl

Concentration: 1 µg/µl

Specificity: Human, mouse: positive

Purity: Affinity purified monoclonal antibody in PBS. Does not contain any preservative.

Storage: Store at -20°C; for long storage, store at -80°C.

Avoid multiple freeze-thaw cycles.

Precautions: This product is for research use only.

Not for use in diagnostic or therapeutic procedures

Description: Monoclonal antibody raised in mouse against Huntingtin, using a recombinant protein.

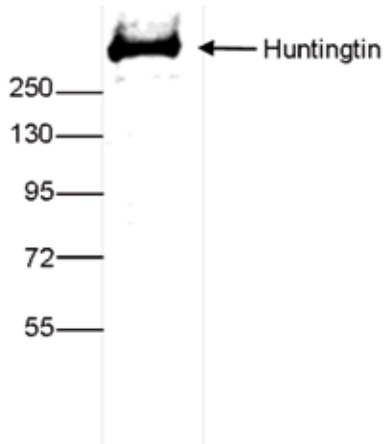
Applications

Applications	Suggested dilution/amount	References
Western blotting	1:1,000	Fig 1
IF	1:400	Fig 2

Target description

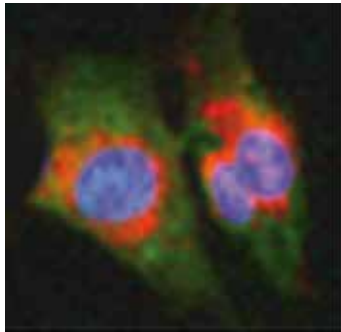
Huntingtin (HTT, UniProt/Swiss-Prot entry P42858) may play a role in microtubule-mediated transport or vesicle function and is required for normal development. It is widely expressed as 2 alternatively polyadenylated transcripts, the larger one being predominantly present in adult and fetal brain. The HTT gene is linked to Huntington's disease, a neurodegenerative disorder characterized by loss of striatal neurons, which is thought to be caused by an expanded, unstable trinucleotide repeat. Whether this leads to loss of function or gain of function is not clear. The expanded repeat may increase the affinity of the HTT protein for the huntingtin-associated protein-1, which is highly expressed in brain and seems to inhibit HTT expression.

Results



Western blot analysis using the Diagenode monoclonal antibody directed against Huntingtin

Mouse brain lysates were analysed by Western blot using the Diagenode antibody against Huntingtin (Cat. No. C15200226) diluted 1:1,000. The position of the protein of interest is indicated on the right; the marker (in kDa) is shown on the left



Immunofluorescence using the Diagenode monoclonal antibody directed against Huntingtin

SKNSH cells were stained with the Diagenode antibody against Huntingtin (Cat. No. C15200226) diluted 1:400 followed by an anti-mouse antibody conjugated to DyLight 488 (green color). WGA conjugated to Alexa Fluor 594 was used to stain the plasma membranes (red). The nuclei were stained with DAPI (blue).

Diagenode sa. BELGIUM | EUROPE

LIEGE SCIENCE PARK
Rue Bois Saint-Jean, 3
4102 Seraing (Ougrée) - Belgium
Tel: +32 4 364 20 50
Fax: +32 4 364 20 51
orders@diagenode.com
info@diagenode.com

Diagenode Inc. USA | NORTH AMERICA

400 Morris Avenue, Suite 101
Denville, NJ 07834 - USA
Tel: +1 862 209-4680
Fax: +1 862 209-4681
orders.na@diagenode.com
info.na@diagenode.com

Last update: September 17, 2015