

DDB2 monoclonal antibody

Other names: DDBB, UV-DDB2

Cat. No. C15200222

Type: Monoclonal

Source: Mouse

Lot #: 001

Size: 50 µg/ 50 µl

Concentration: 1 µg/µl

Specificity: Human: positive

Other species: not tested

Purity: Protein A purified monoclonal antibody in PBS containing 0.05% azide.

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 human DDB2 (Damage-Specific DNA Binding Protein 2) using a KLH-conjugated synthetic peptide containing a sequence from the N-terminal part of the protein.

Applications

	Suggested dilution	Results
Western blotting	1:1,000 - 1:2,000	Fig 1

Target description

DDB2 (UniProt/Swiss-Prot entry Q92466) together with DDB1 forms the heterodimeric DNA damage-binding (DDB) complex. This protein complex binds to DNA following UV damage and functions in nucleotide-excision repair. The DDB complex probably recognizes UV-induced DNA damage and recruits proteins of the nucleotide excision repair (NER) pathway to initiate DNA repair. DDB also mediates the ubiquitylation of histones H3 and H4, which facilitates the cellular response to DNA damage. Defective activity of the DDB complex causes XPE (xeroderma pigmentosum complementation group E), an autosomal recessive disorder characterized by photosensitivity and early onset of carcinomas.

Results

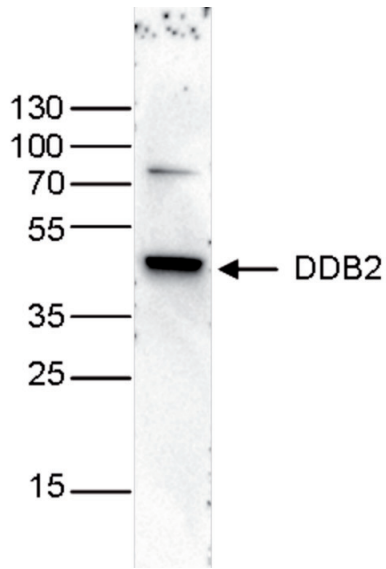


Figure 1. Western blot analysis using the Diagenode monoclonal antibody directed against DDB2

Whole cell extracts from human HeLa cells (25 µg) were analysed by Western blot using the Diagenode antibody against DDB2 (Cat. No. C15200222) diluted 1:1,000 in PBS containing 10% milk. The position of the protein of interest (expected MW 48 kDa) is indicated on the right; the marker (in kDa) is shown on the left.

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: January 15, 2015