

DDB1 antibody (pAb)

Catalog Nos: 61127, 61128

RRID: AB_2793518

Isotype: IgG

Application(s): WB

Reactivity: Human

Volumes: 100 µl, 10 µl

Purification: Affinity Purified

Host: Rabbit

Molecular Weight: 140 kDa

Background: DDB1 (DNA damage-binding protein 1, HBV X-associated protein 1, XAP-1, Xeroderma pigmentosum group E-complementing protein) is a DNA binding protein that associates with DNA following UV damage and functions in excision repair. In addition to its function in DNA repair, DDB1 plays a role in normal mitotic exit by regulating Anaphase Promoting Complex (APC) activity. Depletion of DDB1 in mitotic cells significantly delays mitotic exit. DDB1 facilitates the binding of substrates to the Cul4-Roc1 complex and the ubiquitination of selected proteins, including histones H2A, H3 and H4 localized to the sites of DNA damage. It is also involved in PCNA-dependent polyubiquitination of CDT1 and MDM2-dependent ubiquitination of P53 in response to radiation-induced DNA damage and during DNA replication. In mouse studies, it was identified that loss of DDB1 function prevents hepatocytes from replicating DNA, inducing compensatory proliferation of normal hepatocytes that leads to development of hepatocellular carcinoma.

Immunogen: This DDB1 antibody was raised against a peptide from the C-terminus of human DDB1.

Buffer: Purified IgG in PBS with 30% glycerol and 0.035% sodium azide. Sodium azide is highly toxic.

Application Notes:

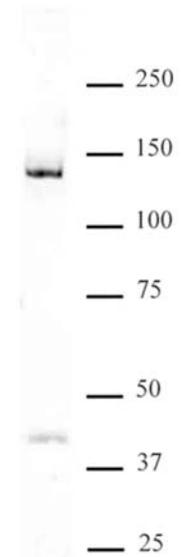
Applications Validated by Active Motif:

WB*: 1:500 - 1:2,000 dilution

*Note: many chromatin-bound proteins are not soluble in a low salt nuclear extract and fractionate to the pellet. Therefore, we recommend a High Salt / Sonication Protocol when preparing nuclear extracts for Western blot.

Storage and Guarantee: Some products may be shipped at room temperature. This will not affect their stability or performance. Avoid repeated freeze/thaw cycles by aliquoting items into single-use fractions for storage at -20°C for up to 2 years. Keep all reagents on ice when not in storage. This product is guaranteed for 12 months from date of receipt.

This product is for research use only and is not for use in diagnostic procedures.



DDB1 antibody (pAb) tested by Western blot
Nuclear extract of HeLa cells (30 µg) probed with DDB1 antibody (1:500).