

D4-GDI antibody (mAb)

Catalog No: 40941

RRID: AB_2793453

Clone: 97A1015

Application(s): WB

Reactivity: Human

Quantity: 100 µg

Purification: Affinity Purified

Host: Mouse

Isotype: IgG1

Concentration: 0.5 µg/µl

Molecular Weight: 23 kDa

Background: D4-GDI (Rho GDP dissociation inhibitor beta or ARHGDI) plays an important role in the activation of the superoxide (O₂⁻)-generating NADPH oxidase of phagocytes. The mechanisms involving cleavage of D4-GDI (Rho GDP) with apoptosis are not presently known. However, activation of the Jun N-terminal kinase, a regulator of apoptosis, may be one of the mechanisms.

Immunogen: This D4-GDI antibody was raised against a synthetic peptide corresponding to amino acid residues in the Fas-induced cleavage site of human D4-GDI.

Buffer: PBS containing 0.02% sodium azide. Sodium azide is highly toxic.

Application Notes:

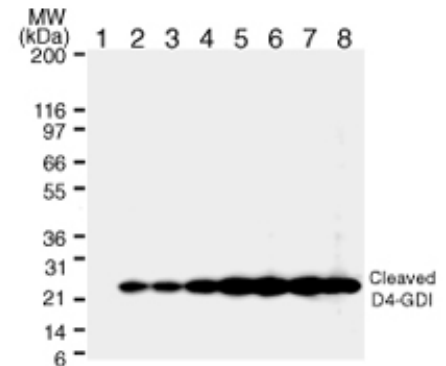
Applications Validated by Active Motif:

WB: 1 - 2 µg/ml dilution

For optimal results, primary antibody incubations should be performed at room temperature. The addition of 0.1% Tween 20 to all blocking solutions may also reduce background. Individual optimization may be required.

Storage and Guarantee: Some products may be shipped at room temperature. This will not affect their stability or performance. Store at 4°C for short term. 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.



D4-GDI mAb tested by Western blot.

Detection of D4-GDI by Western blot analysis. HL-60 cells were treated for different time periods with 10 ng/ml of Fas monoclonal antibody. Harvested cells were lysed, resolved by SDS-PAGE and transferred onto a PVDF membrane. The membrane was incubated with the D4-GDI mAb and immunoreactivity was detected by chemiluminescence. The data shows that the D4-GDI antibody specifically detects the 23 kDa cleaved form of D4-GDI in a time-sensitive manner.