

PRODUCT NAME		
PHF8 polyclonal antibody		
Other names: JHDM1F, MRXSSD, ZNF422		
Cat. No. C15410064 (pAb-064-050)	Type: Polyclonal	Size: 50 µg /39 µl
Lot #: A138-0031	Source: Rabbit	Concentration: 1.3 µg/µl

Description: Polyclonal antibody raised in rabbit against human PHF8 (PHD finger protein 8), using 3 different KLH-conjugated synthetic peptides.

Specificity: Human: positive
Other species: not tested

Applications	Suggested dilution	References
ELISA	1:100 – 1:500	Fig 1
Western blotting	1:1,000	Fig 2

Purity: Affinity purified polyclonal antibody in PBS containing 0.05% azide and 0.05% ProClin 300.

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.

Last data sheet update: March 3, 2010

Target description

PHD finger protein 8 (UniProtKB/Swiss-Prot entry Q9U PP1) is a jumonji domain containing protein. Like other members of the jumonji family, PHF8 may therefore play a role in histone demethylation. Mutations in PHF8 lead to Siderius type X-linked mental retardation (MRXSSD), a mild to borderline type of mental retardation.

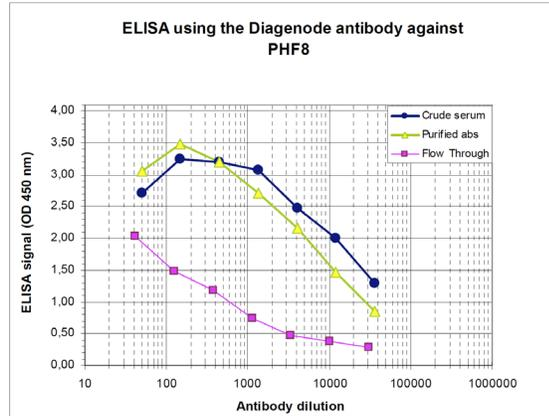


Figure 1
Determination of the antibody titer

To determine the titer of the antibody, an ELISA was performed using a serial dilution of the Diagenode antibody directed against human PHF8 (Cat. No. pAb-064-050), crude serum and Flow Through in antigen coated wells. By plotting the absorbance against the antibody dilution (Figure 1), the titer of the antibody was estimated to be 1:6,500.

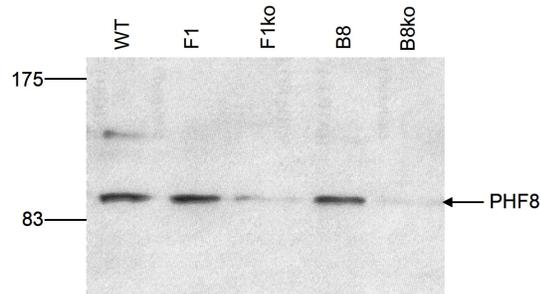


Figure 2
Western blot analysis using the Diagenode antibody directed against PHF8

E14TG2A mouse embryonic stem cells were transfected with a conditional allele of PHF8. Nuclear extracts (135 µg) from wild type cells (WT) and from 2 clones (F1, B8) with an active and a targeted allele (F1ko, B8ko), respectively, were analysed by Western blot using the Diagenode antibody against PHF8 (Cat. No. pAb-064-050). The antibody was 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.

Western blot analysis was performed by Philipp Seibert, BIOTEC, Dept. of Genomics, Prof. A. F. Stewart, TU Dresden, Tatzberg 47/49, 01307 Dresden.