

MORF4 antibody (pAb)

Rabbit Anti-Human/Mouse/Rat MORF4

Instruction Manual

Catalog Number	PK-AB718-4275
Synonyms	MORF4 Antibody: MORF4, Mortality factor 4, cellular senescence-related protein 1, SEN1
Description	Cellular senescence is the terminal non-dividing state that normal cells enter following completion of their proliferative potential. Fusions of immortal human cell lines with each other have led to their assignment to one of four complementation groups. Mortality factor 4 (MORF4) was identified as the lead member of a family of transcription factor-like proteins that reverses this immortal phenotype. Like other members in this family, MORF4 is localized to the nucleus and possesses transcription factor-like motifs such as helix-loop-helix and a leucine zipper motif that might allow it to form transcriptionally active homo- or heterodimers. MORF4 has been shown to bind to the transcription corepressors mSin3A and TLE, suggesting that together, these complexes may play a role in transcriptional repression of genes that lead to cellular senescence.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	MORF4 antibody was raised in rabbits against an 18 amino acid peptide from near the amino terminus of human MORF4.
Purification Method	Affinity chromatography purified via peptide column.
Clone / IgG Subtype	Polyclonal antibody
Species Reactivity	Human, Mouse, Rat
Specificity	
Formulation	Antibody is supplied in PBS containing 0.02% sodium azide.
Reconstitution	During shipment, small volumes of antibody will occasionally become entrapped in the seal of the product vial. For products with volumes of 200 µl or less, we recommend gently tapping the vial on a hard surface or briefly centrifuging the vial in a tabletop centrifuge to dislodge any liquid in the container's cap.
Storage & Stability	Antibody can be stored at 4°C for three months and at -20°C for up to one year. As with all antibodies care should be taken to avoid repeated freeze thaw cycles. Antibodies should not be exposed to prolonged high temperatures.
Applications	E, WB, IHC, IF INote: Antibody might be suitable for other applications not tested so far. Optimal concentrations for each application have to be determined individually. MORF4 antibody can be used for the detection of MORF4 by Western blot at 1 and 2 µg/mL. Antibody can also be used for immunohistochemistry starting at 5 µg/mL. For immunofluorescence start at 20 µg/mL.
Images	Available upon request.
References	Pereira-Smith OM and Smith JR. Genetic analysis of indefinite division in human cells: identification of four complementation groups. Proc. Natl. Acad. Sci. USA 1988; 85:6042-6. Bertram MJ, Berube NG, Hang-Swanson X, et al. Identification of a gene that reverses the immortal phenotype of a subset of cells and is a member of a novel family of transcription factor-like genes. Mol. Cell. Biol. 1999; 1479-85. Yochum GS and Ayer DE. Role for the mortality factors MORF4, MRGX, and MRG15 in transcriptional repression via associations with Pfl1, mSin3A, and transducin-like enhancer of split. Mol. Cell. Biol. 2002; 22:7868-76.
Images	Available upon request.
Related Products	Cat.No. PK-AB718-4275P; MORF4 Peptide Cat.No. PK-AB718-1204; K562 Cell Lysate

FOR IN VITRO RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC PROCEDURES.