

Instruction Manual

Catalog Number	PK-AB718-4031
Synonyms	Emerin Antibody: Emerin, Emery-Dreifuss muscular dystrophy, EDMD, EMD
Description	Emerin is a serine-rich nuclear membrane protein and a member of the nuclear lamina-associated protein family that includes proteins such as LAP2 and MAN1. Each family member, including Emerin, has an ~40 amino acid LEM-domains that binds barrier-to-autointegration (BANF1), a conserved chromatin protein that also serves as a host cell component of retroviral integration complexes, including that of HIV. Emerin is anchored at the inner membrane of the nuclear envelope where it binds to nuclear intermediate filaments that are formed by lamin proteins. Dreifuss-Emery muscular dystrophy is an X-linked inherited degenerative myopathy resulting from mutation in the emerin gene.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	Emerin antibody was raised in rabbits against a 19 amino acid peptide from near the amino terminus of human Emerin.
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. Emerin antibody can be used for detection of Emerin by Western blot at 0.5 - 1 µg/mL. Antibody can also be used for immunohistochemistry starting at 2.5 µg/mL. For immunofluorescence start at 10 µg/mL.
Images	Available upon request.
References	Schirmer EC, Florens L, Guan T, et al. Nuclear membrane proteins with potential disease links found by subtractive proteomics. <i>Science</i> 2003; 301:1380-2. Cai M, Huang Y, Ghirlardo R, et al. Solution structure of the constant region of nuclear envelope protein LAP2 reveals two LEM-domain structures: one binds BAF and the other binds DNA. <i>EMBO J.</i> 2001; 20:4399-407. Chen H and Engelman A. The barrier-to-autointegration protein is a host factor for HIV type 1 integration. <i>Proc. Natl. Acad. Sci. USA</i> 1998; 95:15270-4. Hutchison CJ. Lamins: building blocks or regulators of gene expression? <i>Nat. Rev. Mol. Cell Biol.</i> 2002; 3:848-58.
Images	Available upon request.
Related Products	Cat.No. PK-AB718-4031P; Emerin Peptide Cat.No. PK-AB718-1375; Human Skeletal Muscle Tissue Lysate

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