

BANF1 (CT) antibody (pAb)

Rabbit Anti-Human/Mouse BANF1 (CT)

Instruction Manual

Catalog Number	PK-AB718-4017
Synonyms	BANF1 Antibody: Barrier to autointegration factor 1, BAF, breakpoint cluster region protein 1
Description	Barrier-to-autointegration factor 1 (BANF1) is a conserved chromatin protein that non-specifically binds double-stranded DNA. BANF1 also interacts with a family of nuclear proteins that include LAP2, emerin, and MAN1. It is also a host cell component of retroviral pre-integration complexes (PICs), including that of HIV. BANF1 will bind to p55 Gag (the structural precursor of HIV-1 virions) as well as its cleaved product matrix. In addition to being a host cell component of the PIC, it is thought that BANF1 is also present at low levels in incoming virions, and thus might contribute to the assembly or activity of HIV-1 PICs through direct binding to matrix as well as DNA. Despite its predicted molecular weight, BANF1 often migrates at a higher size in SDS-PAGE.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	BANF1 antibody was raised in rabbits against a 15 amino acid peptide from near the carboxy terminus of human BANF1.
Purification Method	Affinity chromatography purified via peptide column.
Clone / IgG Subtype	Polyclonal antibody
Species Reactivity	Human, Mouse
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, IF Note: Antibody might be suitable for other applications not tested so far. Optimal concentrations for each application have to be determined individually. BANF1 antibody can be used for detection of BANF1 by Western blot at 0.5 - 1 µg/mL. Despite its predicted molecular weight, BANF1 often migrates at a higher size in SDS-PAGE. Antibody can also be used for immunofluorescence starting at 20 µg/mL.
Images	Available upon request.
References	Furukawa K. LAP2 binding protein 1 (L2BP1/BAF) is a candidate mediator of LAP2-chromatin interaction. <i>J. Cell Sci.</i> 1999; 112:2485-92. Cai M, Huang Y, Ghirlando 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.
Images	Available upon request.
Related Products	Cat.No. PK-AB718-4017P; BANF1 Peptide Cat.No. PK-AB718-1305; Human Kidney Tissue Lysate

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