

Bst2 (NT) antibody (pAb)

Rabbit Anti-Human/Mouse Bst2 (Bone marrow stromal cell antigen 2, CD317)

Instruction Manual

Catalog Number	PK-AB718-4661
Synonyms	Bst2 Antibody; Bone marrow stromal cell antigen 2, CD317, HM1.24, tetherin
Description	Bst2 was first identified as a novel terminal B-cell-restricted antigen termed HM1.24 that is expressed on certain bone marrow cells as well as other tissues. The surface expression of Bst2 on fibroblast cell lines facilitated the stromal cell-dependent growth of DW34, a pre-B-cell line, suggesting that this protein is involved in pre-B-cell growth. Later experiments also showed it to be upregulated in several myeloma cells. More recently, Bst2 was identified an inhibitor of retroviral release from human cells whose activity is antagonized by the HIV-1 accessory protein Vpu. While Bst2 caused retention of virions on cell surfaces or endocytosis into Bst2-positive compartments, its depletion abolished the viral requirement for Vpu for virus release. This activity may represent a potential therapeutic strategy for the treatment of HIV/AIDS.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	Nanos3 antibody was raised in rabbits against a 16 amino acid peptide near the carboxy terminus of the human Nanos3.
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 INote: Antibody might be suitable for other applications not tested so far. Optimal concentrations for each application have to be determined individually. Bst2 antibody can be used for detection of Bst2 by Western blot at 1 - 2 µg/mL.
Images	Available upon request.
References	Goto T, Kennel SJ, Abe M, et al. A novel membrane antigen selectively expressed on terminally differentiated human B cells. <i>Blood</i> 1994; 84:1922-30. Ishikawa J, Kaisho T, Tomizawa H, et al. Molecular cloning and chromosomal mapping of a bone marrow stromal cell surface gene, BST2, that may be involved in pre-B-cell growth. <i>Genomics</i> 1995; 26:527-34. Ohtomo T, Sugamata Y, Ozaki Y, et al. Molecular cloning and characterization of a surface antigen preferentially overexpressed on multiple myeloma cells. <i>Biochem. Biophys. Res. Commun.</i> 1999; 258:583-91. Neil SJ, Zang T, and Bieniasz PD. Tetherin inhibits retrovirus release and is antagonized by HIV-1 Vpu. <i>Nature</i> 2008; 451:425-30.
Images	Available upon request.
Related Products	Cat.No. PK-AB718-4661P; Bst2 Peptide Cat. No. PK-AB718-1224; Daudi Cell Lysate

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