

TRIM5 δ antibody (pAb)

Rabbit Anti-Human TRIM5 δ (Retroviral Restriction Factor)

Instruction Manual

Catalog Number	PK-AB718-3241
Synonyms	TRIM5 delta Antibody: TRIM5 delta, TRIM5d
Description	TRIM5 is a member of a broad family of otherwise unrelated proteins defined by the presence of a tripartite motif containing a RING domain, a B-box type 1, and a B-box type 2, followed by a coiled-coil region. TRIM5 has five alternately spliced isoforms, the longest of which is the α -variant which also contains a carboxy-terminal B30.2 (SPRY) domain. Expression of TRIM5 α variants from humans, rhesus monkeys, and African green monkeys enabled resistance to infection by various retroviruses including HIV-1, albeit at differing efficiencies. The TRIM5 δ isoform has been shown to colocalize with the topoisomerase I-interacting proteins BTBD1 and BTBD2 in punctate or elongated cytoplasmic bodies in several mouse and human cells where it appears to serve as a scaffold for the assembly of endogenous BTBD proteins. TRIM5 δ also exhibits autoubiquitination activity in a RING finger- and UbcH5B-dependent manner.
Quantity	100 μ g
Source / Host	Rabbit
Immunogen	Rabbit polyclonal TRIM5 δ antibody was raised against a synthetic peptide corresponding to 13 amino acids near the carboxy-terminus of human TRIM5 δ (Genbank accession No. NP_149084).
Purification Method	Affinity chromatography purified via peptide column.
Clone / IgG Subtype	Polyclonal antibody
Species Reactivity	Human
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 μ l 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.h application have to be determined individually. TRIM5 delta antibody can be used for detection of TRIM5 delta by Western blot at 2 μ g/mL. A band at approximately 25 kDa can be detected. Antibody can also be used for immunohistochemistry starting at 1 μ g/mL. For immunofluorescence start at 10 μ g/mL.
Images	Available upon request.
References	Reymond A, Meroni G, Fantozzi A, et al. The tripartite motif family identifies cell compartments. EMBO J. 2001; 20:2140-51. Stremmlau M, Owens CM, Perron MJ, et al. The cytoplasmic body component TRIM5 δ restricts HIV-1 infection in Old World monkeys. Nature 2004; 427:848-53. Hatziioannou T, Perez-Caballero D, Yang A, et al. Retrovirus resistance factors REF1 and Lv1 are species-specific variants of TRIM5 α . Proc. Nat'l. Acad. Sci. USA 2004; 101:10774-9 Xu L, Yang L, Moitra PK, et al. BTBD1 and BTBD2 colocalize to cytoplasmic bodies with the RBCC/tripartite motif protein, TRIM5 δ . Exp. Cell Res. 2003; 288:84-93.
Images	Available upon request.
Related Products	Blocking Peptide, Cat. No. PK-AB718-3241P; Mouse brain Cell Lysate, Cat. No. PK-AB718-1403 TRIM5-alpha Antibody (IN1), Cat. No. PK-AB718-3247; TRIM5-alpha Antibody (IN2), Cat. No. PK-AB718-3251 TRIM5-alpha Antibody (CT), Cat. No. PK-AB718-3249

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