

# TRIM5 $\gamma$ antibody (pAb)

## Rabbit Anti-Human TRIM5 $\gamma$ (Retroviral Restriction Factor)

### Instruction Manual

|                     |   |
|---------------------|---|
| Catalog Number      | PK-AB718-3239   |
| Synonyms            | TRIM5 gamma Antibody; TRIM5 gamma, TRIM5g   |
| 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 $\alpha$ -variant which also contains a carboxy-terminal B30.2 (SPRY) domain. While a function has not yet been assigned to TRIM5 $\gamma$ , it is known that expression of TRIM5 $\gamma$ variants from humans, rhesus monkeys, and African green monkeys enable resistance to infection by various retroviruses including HIV-1, albeit at differing efficiencies. Furthermore, the TRIM5 $\gamma$ isoform appears to serve as a scaffold for the assembly of endogenous BTBD1 and BTBD2 proteins and also exhibits autoubiquitination activity in a RING finger- and UbCH5B-dependent manner. |
| Quantity            | 100 $\mu$ g   |
| Source / Host       | Rabbit  |
| Immunogen           | Rabbit polyclonal TRIM5 $\gamma$ antibody was raised against a synthetic peptide corresponding to 15 amino acids near the carboxy-terminus of human TRIM5 $\gamma$ (Genbank accession No. NP_149083).   |
| 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 $\mu$ 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 <span style="float: right;">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.</span><br>TRIM5 gamma antibody can be used for detection of TRIM5 gamma by Western blot at 2 $\mu$ g/mL. A band at approximately 35 kDa can be detected. Antibody can also be used for immunohistochemistry starting at 2 $\mu$ g/mL. For immunofluorescence start at 20 $\mu$ 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.<br>Stremlau M, Owens CM, Perron MJ, et al. The cytoplasmic body component TRIM5 $\gamma$ restricts HIV-1 infection in Old World monkeys. Nature 2004; 427:848-53.<br>Hatzioannou T, Perez-Caballero D, Yang A, et al. Retrovirus resistance factors REF1 and Lv1 are species-specific variants of TRIM5 $\alpha$ . Proc. Nat'l. Acad. Sci. USA 2004; 101:10774-9<br>Xu L, Yang L, Moitra PK, et al. BTBD1 and BTBD2 colocalize to cytoplasmic bodies with the RBCC/tripartite motif protein, TRIM5 $\gamma$ . Exp. Cell Res. 2003; 288:84-93.   |
| Images              | Available upon request.   |
| Related Products    | Blocking Peptide, Cat. No. PK-AB718-3239P; Human bladder Cell Lysate, Cat. No. PK-AB718-1310<br>TRIM5-alpha Antibody (IN2), Cat. No. PK-AB718-3251; TRIM5-alpha Antibody, Cat. No. PK-AB718-3241<br>TRIM5-alpha Antibody (IN1), Cat. No. PK-AB718-3247; TRIM5-alpha Antibody (CT), Cat. No. PK-AB718-3249   |

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