

# TRIM25 (CT) antibody (pAb)

Rabbit Anti-Human/Mouse TRIM25 (CT)

## Instruction Manual

Catalog Number	PK-AB718-4315
Synonyms	TRIM25 Antibody: Tripartite motif-containing 25, estrogen responsive finger protein, EFP
Description	TRIM25, also known as estrogen responsive finger protein (EFP) is a member of the RING finger-B box-coiled coil family and is a downstream target of estrogen receptor $\alpha$ . Mice lacking this gene displayed an underdeveloped uterus and reduced estrogen responsiveness demonstrating that this protein is essential for estrogen dependent proliferation. TRIM25 expression is increased in breast carcinoma where it promotes the growth of breast tumor by functioning as an E3 ubiquitin ligase that targets the negative cell cycle checkpoint 14-3-3s. Recent reports have indicated that TRIM25 E3 ubiquitin ligase is essential for the retinoic-acid-inducible gene 1 (RIG-1) signaling pathway to elicit host antiviral innate immunity.
Quantity	100 $\mu$ g
Source / Host	Rabbit
Immunogen	TRIM25 antibody was raised in rabbits against a 19 amino acid peptide from near the carboxy terminus of human TRIM25.
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 $\mu$ 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, ICC, IF <b>INote:</b> Antibody might be suitable for other applications not tested so far. Optimal concentrations for each application have to be determined individually. TRIM25 antibody can be used for detection of TRIM25 by Western blot at 1 $\mu$ g/mL. Antibody can also be used for immunocytochemistry starting at 20 $\mu$ g/mL. For immunofluorescence start at 20 $\mu$ g/mL.
Images	Available upon request.
References	Inoue S, Orimo A, Hosoi T, et al. Genomic binding-site cloning reveals an estrogen-responsive gene that encodes a RING finger protein. Proc. Natl. Acad. Sci. USA 1993; 111:17-20. Orimo A, Inoue S, Minowa O, et al. Underdeveloped uterus and reduced estrogen responsiveness in mice with disruption of the estrogen-responsive finger protein gene, which is a direct target of estrogen receptor $\alpha$ . Proc. Natl. Acad. Sci. USA 1999; 96:12027 Thomson SD, Ali S, Pickles L, et al. Analysis of estrogen-responsive finger protein expression in benign and malignant human breast. Int. J. Cancer 2001; 91:152-8. Urano T, Saito T, Tsukui T, et al. Efp targets 14-3-3s for proteolysis and promotes breast tumour growth. Nature 2002; 417:871-5.
Images	Available upon request.
Related Products	Cat.No. PK-AB718-4315P; TRIM25 Peptide Cat.No. PK-AB718-1201; HeLa Cell Lysate

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