

EVER2 (IN) antibody (pAb)

Rabbit Anti-Human/Mouse/Rat EVER2 (Transmembrane channel-like protein 8, TMC8)

Instruction Manual

Catalog Number	PK-AB718-4605
Synonyms	EVER2 Antibody: Transmembrane channel-like protein 8, TMC8
Description	Epidermodysplasia verruciformis (EV) is an autosomal recessive genodermatosis associated with a high risk of skin cancers resulting from a high susceptibility to infection by specific human papillomaviruses. Mutations in two homologous genes EVER1 and EVER2 cause the majority of EV cases. These two proteins form a complex and interact with the zinc transporter ZnT-1 in the endoplasmic reticulum. Cells lacking EVER2 accumulated higher levels of zinc in the nucleolus and nucleus compare to those cells with and intact EVER2 gene, indicating that one role of EVER2 is to regulate the intracellular distribution of zinc. At least three isoforms of EVER2 are known to exist; this antibody will only recognize the larger isoform. EVER2 has no cross-reactivity to EVER1.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	EVER2 antibody was raised in rabbits against a 14 amino acid peptide from near the center of human EVER2.
Purification Method	Affinity chromatography purified via peptide column.
Clone / IgG Subtype	Polyclonal antibody
Species Reactivity	Human, Mouse, Rat
Specificity	At least three isoforms of EVER2 are known to exist; this antibody will only recognize the larger isoform. EVER2 has no cross-reactivity to EVER1.
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, IHC, IF INote: Antibody might be suitable for other applications not tested so far. Optimal concentrations for each application have to be determined individually. EVER2 antibody can be used for detection of EVER2 by Western blot at 1 - 2 µg/mL. Antibody can also be used for immunohistochemistry starting at 5 µg/mL. For immunofluorescence start at 20 µg/mL.
Images	Available upon request.
References	Ramoz N, Taieb A, Rueda L-A, et al. Evidence for a nonallelic heterogeneity of epidermodysplasia verruciformis with two susceptibility loci mapped to chromosome regions 2p21-p24 and 17q25. <i>J. Invest. Dermatol.</i> 2000; 114:1148-53. Ramos N, Rueda L-A, Bouadjar B, et al. Mutations in two adjacent novel genes are associated with epidermodysplasia verruciformis. <i>Nat. Genet.</i> 2002; 32:579-81. Lazarczyk M, Pons C, Mendoza J-A, et al. Regulation of cellular zinc balance as a potential mechanism of EVER-mediated protection against pathogenesis by cutaneous oncogenic human papillomaviruses. <i>J. Exp. Med.</i> 2008; 205:35-42.
Images	Available upon request.
Related Products	Cat.No. PK-AB718-4605P; EVER2 Peptide Cat. No. PK-AB718-1205; Jurkat Cell Lysate

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