

Instruction Manual

Catalog Number	PK-AB718-4053
Synonyms	VISA Antibody: Virus-induced signaling adapter, mitochondrial antiviral signaling protein, MAVS, CARD adapter inducing interferon-beta, Cardif, IPS-1
Description	Two distinct signaling pathways activate the host innate immunity against viral infection. One pathway is reliant on members of the Toll-like receptor (TLR) family while the other uses the RNA helicase RIG-I as a receptor for intracellular viral double-stranded RNA as a trigger for the immune response. VISA is a mitochondrial membrane protein that was identified as a critical component in the IFN- β signaling pathways that recruits IRF-3 to RIG-I, leading to its activation and that of NF- κ B. VISA is also thought to interact with other components of the innate immune pathway such as the TLR adapter protein TRIF, TRAF2 and TRAF6. VISA also interacts with the IKK α , IKK β and IKK ϵ kinases through its C-terminal region. Cleavage of this region by the Hepatitis C virus (HCV) protease allows HCV to escape the host immune system. At least three isoforms of VISA are known to exist.
Quantity	100 μ g
Source / Host	Rabbit
Immunogen	VISA antibody was raised in rabbits against a 13 amino acid peptide from near the amino terminus of human VISA.
Purification Method	Affinity chromatography purified via peptide column.
Clone / IgG Subtype	Polyclonal antibody
Species Reactivity	Human, Mouse, Rat
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. VISA antibody can be used for detection of VISA by Western blot at 0.5 - 2 μ g/mL. Antibody can also be used for immunohistochemistry starting at 5 μ g/mL. For immunofluorescence start at 20 μ g/mL.
Images	Available upon request.
References	Seth RB, Sun L, and Chen ZJ. Antiviral innate immunity pathways. Cell Res. 2006; 16:141-7. Xu LG, Wang YY, Han KJ, et al. VISA is an adapter protein required for virus-triggered IFN- β signaling. Mol. Cell 2005; 19:727-40. Meylan E, Curran J, Hofman K, et al. Cardif is an adaptor protein in the RIG-I antiviral pathway and is targeted by hepatitis C virus. Nature 2005; 1167-72.
Images	Available upon request.
Related Products	Cat.No. PK-AB718-4053P; VISA Peptide Cat.No. PK-AB718-1288; A20 Cell Lysate

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