

HCAM Polyclonal antibody (pAb)

Rabbit Anti-Human/Mouse/Rat Homing Cell Adhesion Molecule (CD44)

Instruction Manual

Catalog Number	PK-AB577-3393
Description	Cell adhesion molecules (CAMs) are a family of closely related cell surface glycoproteins involved in cell-cell interactions during growth and are thought to play an important role in embryogenesis and development. The CAM family includes NCAM (neuronal cell adhesion molecule), ICAM (intracellular adhesion molecule), PECAM (platelet/endothelial cell adhesion molecule), BLCAM (produced by alternative splicing of the CD22 gene), VCAM (vascular cell adhesion molecule), and HCAM (homing cell adhesion molecule, also designed CD44).
Quantity	100 µg
Source / Host	Rabbit
Immunogen	Synthetic peptide surrounding amino acid 39 of human HCAM/CD44 (ID BV-376)
Clone / IgG Subtype	Polyclonal antibody
Species Reactivity	Human, Mouse, Rat
Specificity	NA
Formulation	100 µg (0.2 mg/ml) protein A purified rabbit polyclonal antibody in phosphate-buffered saline (PBS) containing 50% glycerol, 1% BSA, and 0.02% thimerosal.
Reconstitution	Can be diluted in other aqueous buffers at the concentrations determined for the respective application.
Storage & Stability	Store at -20°C. For long-term storage, aliquot and freeze at -70°C. Avoid repeated freeze/thaw cycles.
Applications	The antibody has been used in Western blotting (0.5-4 µg/ml). Based on customer's feedback, the antibody also works well in immunoprecipitation (20 µg/ml) and immunohistochemistry (20 µg/ml). However, the optimal concentrations should be determined individually. Blocking peptide is available separately (Cat.# PK-AB577-3393P).

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