

## Instruction Manual

Catalog Number	PK-AB704-68505
Quantity	100 µg
Clone	H572
IgG Subtype	mouse IgG1
Purity	Protein A affinity chromatography
Specificity	Reacts the glycoprotein EndoGlyx-1 that is expressed on endothelial cells of human normal and tumor blood vessel, but not on cells of mesenchymal, neuroectodermal and epithelial derivation.
Cross-reactivity	There are no data about cross-reactivity with other species.
Formulation	PBS
Storage	Stable at 2-8°C. After opening, aliquote contents and freeze at -20°C for long-term storage. Avoid repeated freeze/thaw cycles.
Shipping conditions	2-8°C
Applications	Flow cytometry Immunoprecipitation Immunohistology:PK-AB704-68505 is suitable for stainings on unfixed, acetone-fixed or ethanol-fixed frozen tissues, but not on formalin-fixed, paraffin-embedded archival material. Acetone-fixed frozen tissues showed optimal staining results. PK-AB704-68505 is not suitable for Western Blot analysis. Optimal antibody concentrations should be determined by the individual laboratory for each application.

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