

Instruction Manual

Catalog Number	PK-AB704-68530
Quantity	100 µg
Clone	F11-24
IgG Subtype	mouse IgG1
Purity	Protein A affinity chromatography
Specificity	Reacts with Fibroblast Activation Protein (FAP), a cell surface glycoprotein of reactive human fibroblast. Among the normal adult organs tested PK-AB704-68530 does not show immunoreactivity in immunohistological studies with either resting fibroblasts (fibrocytes) of the normal organ mesenchyme or the parenchymal cells. The single exception noted is a subset of endocrine cells in the pancreatic islets, presumably glucagon-producing A cells.
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-AB704-68530 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. 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.