

Instruction Manual

Catalog Number	PK-AB714-64050HM
Quantity	100 µg
Description	Monoclonal antibodies were produced by immunizing Syrian hamsters with a murine thymic stromal cell line. This antibody recognizes mouse Podoplanin. Hamster IgG antibody from hybridomas was purified from cell culture supernatant by Protein G chromatography.
Host species	Syrian Hamster
Antigen	NA
Stabilizer	none
Purification	Protein G affinity chromatography
Specificity	The monoclonal antibody will detect Podoplanin on the surface of mouse lymphatic endothelial cells and some epithelial cell types by immunostaining or immunohistochemistry. It works well with paraffin sections. The antibody will detect a 36-40 kDa band in Western blotting. It also works with immunoprecipitation:
Formulation	PBS (pH 7.4) w/o preservative
Reconstitution	When reconstituted in sterile water to a concentration of 1.0 mg/ml the antibody is stable for at least six weeks at 2-4°C.
Storage & Stability	The lyophilized antibody, thought stable at room temperature, is best stored desiccated below 0°C. Reconstituted antibody is stable at 4°C for > one month or can be stored in working aliquots at -20°C for more than six months. Cross-Reactivity: The antibody does not cross-react with human Podoplanin. Other species not tested.
Applications	Immunostaining/Immunohistochemistry: Use diluted 1-10 µg/ml IgG. Western blotting: Use 1-10 µg/ml IgG. Immunoprecipitation: For IP use 1-5 µg IgG per 1 ml lysate or reaction volume.
References	Schacht et al., EMBO J 22:3546, 2003

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