

# Podoplanin antibody (mAb)

PromoKine

Mouse Anti-Human Podoplanin (Clone 4D5aE5E6) TNF Family

## Instruction Manual

Catalog Number	PK-AB714-64050
Quantity	100 µg
Description	Monoclonals were produced with the help of BALB/c mice using a membrane protein fraction of isolated human glomeruli as the immunizing antigen. The mouse IgG1kappa antibody (#4D5aE5E6) from hybridomas was purified from cell culture supernatant by Protein G chromatography.
Host species	Mouse
Antigen	Membrane protein fraction of isolated human glomeruli
Stabilizer	none
Purification	Protein G affinity chromatography
Specificity	The antibody is human-specific.
Formulation	Lyophilized from solution in PBS (pH 6.0) w/o preservative
Reconstitution	The lyophilized IgG is stable at 4°C. for at least one month and for more than a year when kept at -20°C. When reconstituted in sterile water to a concentration of 0.1-1.0 mg/ml the antibody is stable for at least six weeks at 2-4°C in the presence of a preservative. Avoid repeated freeze-thaw cycles.
Storage & Stability	Reconstituted antibody can also be aliquotted and stored at -20°C to -70°C for at least 6 months without detectable loss of activity.
Applications	Immunohistochemistry/Immunofluorescence: Cryo sections (acetone fixed); Paraffin sections (after autoclave or microwave treatment in 0.01 M citrate buffer, pH 6.0). Western blotting: Use at 1-5 µg/ml IF/IHC: Use at 1-2 µg/ml Optimal dilutions should be determined by each laboratory for each application.
References	Breiteneder-Geleff et al., Am. J. Pathol. 151:1141,1997 Breiteneder-Geleff et al., Am. J. Pathol. 154:385,1999 Kerjaschki et al., Nature Med. 12:230, 2006 Kriehuber et al., J. Exp. Med. 194:797, 2001

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