

Instruction Manual

Catalog Number	PK-AB714-64050MR
Quantity	100 µg
Description	Monoclonals were produced with the help of BALB/c mice using a membrane protein fraction of isolated rat glomeruli as the immunizing antigen. The mouse IgG antibody (#LF3(B7)D5B3) from hybridomas was purified from cell culture supernatant by Protein G chromatography.
Host species	Mouse
Antigen	Membrane protein fraction of isolated rat glomeruli
Stabilizer	none
Purification	Protein G affinity chromatography
Specificity	The antibody is rat specific!
Formulation	PBS w/o preservative
Reconstitution	Centrifuge vial prior to opening. Reconstitute in sterile, ultrapure water to a concentration of 0.1-1.0 mg/ml.
Storage & Stability	The lyophilized antibody is stable at RT for at least one month and for more than a year when kept at -20°C. When reconstituted in sterile water to a concentration, the antibody is stable for at least two weeks at 2-4°C in the presence of a preservative. Frozen aliquots are stable for at least six months at -20°C. Avoid repeated freeze-thaw cycles.
Applications	Immunohistochemistry: Cryo section (acetone-fixed) Western blotting: Use at 1-5 µg/ml. Optimal dilutions should be determined by each laboratory for each application.
References	Breiteneder-Geleff et al., Am. J. Pathol. 151:1141,1997 Schacht et al., EMBO J. 22:3546-56, 2003

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