

Instruction Manual

Catalog Number	PK-AB704-69098F
Quantity	100 tests
Clone	MnS-1
IgG Subtype	mouse IgG1
Purity	> 95% IgG1 as determined by SDS-gel electrophoresis prior to addition of stabilizer and preservative.
Specificity	Specifically recognizes human MnSOD.
Cross-reactivity	There are no data about cross-reactivity with other species.
Formulation	50 mM TRIS 100 mM NaCl 1% BSA 0.02% NaN , pH 7.4
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	Immunohistochemistry: PK-AB704-69098 can be used to stain both indirectly or directly acetone-fixed cryostat sections or cell smears. Recommended for PK-AB704-69098 is the alkaline phosphatase-anti-alkaline-phosphatase (APAAP), or peroxidase anti-peroxidase (PAP) procedures or the three stage immunoperoxidase technique on acetone-fixed cryostat sections. PK-AB704-69098 may be used at a concentration of 0,5-10 µg/ml. Western blot: PK-AB704-69098 is suitable for Western blot application. Under reducing conditions/SDS electrophoresis using BMS122 a 25 kDa band will be recognized (monomer of MnSOD protein). The native form of the protein with a molecular weight of 80 kDa will be recognized under non-reducing conditions. 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.