

Bax antibody (mAb)

Mouse Anti-Human/Mouse/Rat Bax (Bcl-2 Associated X-Protein)

Instruction Manual

Catalog Number	PK-AB704-03352
Quantity	100 µg
Clone	6A7
IgG Subtype	mouse IgG1
Purity	NA
Specificity	Binds to Bax of many different species. 6A7 avidly binds the monomeric form of Bax, but not Bax complexed with either Bcl-XL or Bcl-2.
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	Western Blot: recommended concentration 0.5 - 1 µg/ml ELISA: recommended concentration 0.5 - 1 µg/ml Immunofluorescence: recommended concentration 0.5 - 1 µg/ml Immunoprecipitation in detergent 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.