

# Bax antibody (mAb)

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

### Instruction Manual

Catalog Number	PK-AB704-03351
Quantity	100 µg
Clone	2D2
IgG Subtype	mouse IgG1
Purity	NA
Specificity	Reacts with human Bax.
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 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.