

Instruction Manual

Catalog Number	PK-AB704-69056
Quantity	100 µg
Clone	8C8
IgG Subtype	mouse IgG1
Purity	NA
Specificity	Reacts with human 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. Further applications have not been evaluated, but suitability is expected. 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.