

Instruction Manual

Catalog Number	PK-AB704-69227
Quantity	100 µg
Clone	14C2
IgG Subtype	mouse IgG1
Purity	> 95%
Specificity	Specifically recognizes human Fas Ligand.
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	ELISA Optimal antibody concentrations have to be determined by the individual laboratory for each application.

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