

Bcl-6 antibody (pAb)

Rabbit Anti-Human/Mouse/Rat Bcl-6

Instruction Manual

Catalog Number	PK-AB577-3408
Description	Bcl-6 encodes a protein of 706 amino acids with a predicted molecular weight of 79 kDa. The Bcl-6 protein has homology with zinc finger transcription factors. However, little is known about its functions.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	Synthetic peptide mapping to amino acid residues near the C-terminus of human Bcl-6.
Clone / IgG Subtype	Rabbit IgG
Species Reactivity	human, mouse, rat
Specificity	See Applications or Species Reactivity.
Formulation	100 µg (0.2 mg/ml) affinity purified rabbit anti-Bcl-6 polyclonal antibody in phosphate buffered saline (PBS), pH 7.2, containing 30% glycerol, 0.5% BSA, 0.01% thimerosal.
Reconstitution	Can be diluted in other aqueous buffers at the concentrations determined for the respective application.
Storage & Stability	Store at -20°C. For long-term storage, aliquot and refreeze at -70°C. Avoid repeated freeze/thaw cycles.
Applications	Western blot (0.5-4 µg/ml), Neutralization (0.5-5 µg/ml), immunoprecipitation (10-20 µg/ml), and immunohistochemistry (20-40 µg/ml). The optimal conditions should be determined individually. Antibody might be suitable for other applications not tested so far.

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