

Instruction Manual

Catalog Number	PK-AB577-3805
Description	Prohibitin is an evolutionarily conserved protein that has anti-proliferative activity. Prohibitin is a 30 kDa, post-synthetically modified protein found in the inner membrane of mitochondria and on the plasma membrane of B cells. It mediates B cell mutation. Prohibitin Controls cell cycle in both transcriptional and post-translational mechanisms. Prohibitin is a potential tumor suppressor. Prohibitin gene mutations are correlated with the breast cancer development and/or progression in over 80% of the cell lines analyzed.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	Synthetic peptide surrounding amino acid 256 of human Prohibitin.
Clone / IgG Subtype	Rabbit IgG
Species Reactivity	human, mouse, rat
Specificity	See Applications or Species Reactivity.
Formulation	100 µg (0.5 mg/ml) affinity purified rabbit 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 blotting (0.5-4 µg/ml), immunoprecipitation (10-20 µg/ml) and immunofluorescence (5-10 µg/ml). The optimal conditions should be determined individually. Antibody might be suitable for other applications not tested so far. The antibody detects a 33 kDa protein, corresponding to the apparent molecular mass of Prohibitin on SDS-PAGE immunoblots, in samples from human, mouse and rat origins.

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