

Instruction Manual

Catalog Number	PK-AB577-3773
Description	Steroid receptor coactivator 3, NCOA3, Nuclear Receptor Coactivator 3; TRAM1, Thyroid Hormone Receptor Activator Molecule 1; RAC3, Receptor-Associated Coactivator 3; AIB1, Amplified in Breast Cancer-1; pCIP, CBP-Interactin protein
Quantity	100 µl
Source / Host	Rabbit
Immunogen	Synthetic peptide of human SRC3
Clone / IgG Subtype	Rabbit IgG
Species Reactivity	human
Specificity	See Applications or Species Reactivity.
Formulation	100 µl affinity purified rabbit polyclonal antibody in phosphate-buffered saline (PBS) containing 30% glycerol, 1% BSA and 0.02% 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 (1:500 – 1:2500) and immunoprecipitation. The optimal concentrations should be determined individually. Antibody might be suitable for other applications not tested so far. HeLa cell lysate can be used as a positive control. The antibody recognizes the SRC3 of human origin. Reactivity to other species has not been tested.

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