

# 4-1BBR antibody (pAb)

## Rabbit Anti-Human 4-1BB Receptor

### Instruction Manual

Catalog Number	PK-AB577-5370
Description	Human 4-1BB Receptor (4-1BBR) is an inhibitor of 4-1BBL and can block 4-1BBL bioactivity. Human 4-1BBR is a 17.7 kDa protein containing 167 amino acid residues.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	Recombinant human 4-1BBR (Internal ID BV-808)
Clone / IgG Subtype	Polyclonal antibody
Species Reactivity	Human (Other species have not been tested)
Specificity	NA
Formulation	100 µg (0.5 mg/ml) affinity purified rabbit anti-human 4-1BBR polyclonal antibody in 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 freeze at -70°C. Avoid repeated freezing/defrost cycles.
Applications	The antibody can be used in Western Blot analysis (0.5-2.0 µg/ml). Per researcher's feedback, it can also be used in Neutralization (2-4 µg/ml). However, the optimal conditions should be determined individually.

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