

### Instruction Manual

Catalog Number	PK-AB577-3851
Description	Xenopus Dickkopf (Dkk)-1 was initially discovered as a Wnt antagonist that plays an important role in head formation. By far, four members of Dkk have been identified in mammals. Each Dkk molecule contains two conserved cysteine-rich domains. Recent studies showed that the second Cys-rich domains of Dkk1 and Dkk2 inhibited Wnt-3a-activated signaling, whereas the first Cys-rich domains had no effects. In addition, the second Cys-rich domain of Dkk-2, but not that of Dkk-1, was able to activate the canonical pathway in the presence of LRP6, and this LRP-dependent signaling does not require Dvl.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	Synthetic peptide surrounding amino acid 255 of mouse Dkk1.
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). The optimal conditions should be determined individually. Antibody might be suitable for other applications not tested so far. The antibody detects ~35 kDa Dkk1 of human, mouse and rat origins.

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