

### Instruction Manual

<b>Catalog Number</b>	PK-AB577-3429
<b>Description</b>	DNA mismatch repair genes have been found to be associated with hereditary nonpolyposis colorectal cancer (HNPCC). Inherited mutations in the MSH2 were demonstrated at high frequency in HNPCC and were shown to be associated with microsatellite instability. The demonstration that 10 to 45% of pancreatic, gastric, breast, ovarian and small cell lung cancers also display microsatellite instability suggests that DNA mismatch repair is not restricted to HNPCC tumors but is a common feature in tumor initiation or progression.
<b>Quantity</b>	100 µg
<b>Source / Host</b>	Rabbit
<b>Immunogen</b>	Synthetic peptide mapping to a region between 1-50 of MSH2.
<b>Clone / IgG Subtype</b>	Rabbit IgG
<b>Species Reactivity</b>	human, mouse, rat
<b>Specificity</b>	See Applications or Species Reactivity.
<b>Formulation</b>	100 µg (0.2 mg/ml) purified rabbit polyclonal antibody in phosphate-buffered saline (PBS) containing 50% glycerol, 1% BSA, and 0.02% sodium azide.
<b>Reconstitution</b>	Can be diluted in other aqueous buffers at the concentrations determined for the respective application.
<b>Storage &amp; Stability</b>	Store at -20°C. For long-term storage, aliquot and refreeze at -70°C. Avoid repeated freeze/thaw cycles.
<b>Applications</b>	Western blotting (1-4 µg/ml), immunoprecipitation (15-25 µg/ml), and immunohistochemistry (15-25 µg/ml). The optimal conditions should be determined individually. Antibody might be suitable for other applications not tested so far. The antibody recognizes the MSH2 of human, mouse, and rat origins. Reactivity to other species has not been tested.

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