

**6-His Monoclonal Antibody, Purified**  
**Catalog Number: MMS-156P**  
**Available Size: 0.2 mL, 0.5 mL**

**Description:** Monoclonal Antibody, 6-His (6His, 6xHis, hexaHis)

**Clone:** 6-His

**Form:** Purified Antibody (Stored in PBS + 0.03% Thimerosal)

**Host:** Mouse

**IsoType:** IgG1

**[Ab]:** 1 mg/mL

**Specificity:** Monoclonal 6-His antibody was raised against a recombinant protein containing the sequence HHHHHH. The sequence of the protein flanking the poly His tag does not influence antibody binding.

**Uses:** This antibody is effective in immunoblotting, immunoprecipitation and immunofluorescence.

**Suggested Working Dilution:** The optimal working dilution should be determined for each specific assay condition. *This antibody is sold for laboratory research use only, not for human or in-vivo use. Covance antibodies may not be resold or modified for resale without prior written approval.*

- Western blot: 1:1,000\*
- Immunofluorescence: 1:1,000
- Immunoprecipitation: 1:150

**Notes:** \*Our Posi-Tag Control Protein (PRO-133P) can be used as a helpful positive control

for this antibody.

**Storage:** Store at -20°C. It is suggested that the total volume be divided into useable aliquots upon initial thaw.

**References:** Deinert K, Fasiolo F, Hurt EC, Simos G. Arc1p organizes the yeast aminoacyl-tRNA synthetase complex and stabilizes its interaction with the cognate tRNAs. *J. Biol. Chem.* 276(8):6000-6008, 2001.

Kojima H, Nemoto A, Uemura T, Honma R, Ogura M, Liu Y. rDrak1, a novel kinase related to apoptosis, is strongly expressed in active osteoclasts and induces apoptosis. *J. Biol. Chem.* 276(22):19238-19243, 2001.

Lapinski PE, Miller GG, Tampe R, Raghavan M. Pairing of the nucleotide binding domains of the transporter associated with antigen processing. *J. Biol. Chem.* 275(10):6831-6840, 2000.

Bartsch D, Casadio A, Karl KA, Serodio P, Kandel ER. CREB1 encodes a nuclear activator, a repressor, and a cytoplasmic modulator that form a regulatory unit critical for long-term facilitation. *Cell* 95:211-223, 1998.

**Warranty/Conditions:** Covance products may not be resold or modified for resale without prior written approval.

**Related Products:**

- 6-His Polyclonal Antibody  
Catalog Number PRB-156C
- Posi-Tag Epitope Tag Control Protein  
Catalog Number PRO-133P

- 6-His Monoclonal Antibody  
Catalog Number MMS-156R
  
- 6-His Monoclonal Antibody, Affinity  
Matrix  
Catalog Number AFC-156P

**Product Revision Date: 2/18/2007**