

AU5 Monoclonal Antibody, Purified
Catalog Number: MMS-135P
Available Size: 0.2 mL, 0.5 mL

Description: Monoclonal Antibody, AU5

Intended Use: **Research Use Only (RUO)**

This product is sold for laboratory research use only, not for human or in-vivo use.

Clone: AU5

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

Host: Mouse

IsoType: IgG1

[Ab]: 1 mg/mL

Specificity: The AU5 monoclonal antibody was raised against an SDS-disrupted BPV-1 preparation and mapped to the six amino acid epitope TDFYLK. This clone was specifically designed for use in epitope tagging. Because of its small size, clone AU5 is unlikely to alter the activity of the cloned sequence. This antibody was purified using protein-G chromatography.

Uses: This antibody is effective in immunofluorescence (IF), immunoblotting, and immunoprecipitation (IP) of tagged proteins.

Suggested Working Dilution: The optimal working dilution should be determined for each specific assay condition.

- Western blot: 1:500*
- IF: 1:500

- IP: 1:150

Storage: Store at -20°C. Upon initial thawing, apportion into working aliquots and store at -20°C. Avoid repeated freeze-thaw cycles to prevent denaturing the antibody. For long-term storage, keep the antibody at -80°C.

References: Burke TW, Cook JG, Asano M, Nevins JR. Replication factors MCM2 and ORC1 interact with the histone acetyltransferase HBO1. *J Biol Chem* 276(18):15397-15408, 2001.

Marinissen MJ, Chiariello M, Gutkind JS. Regulation of gene expression by the small GTPase Rho through the ERK6 (p38 gamma) MAP kinase pathway. *Genes Dev* 15(5):535-553, 2001.

Lim PS, Jensen AB, Cowser L, Nakai Y, Lim LY, Jin XW, Sundberg JP. Distribution and specific identification of papillomavirus major capsid protein epitopes by immunocytochemistry and epitope scanning of synthetic peptides. *588-588fect Dis* 162-1269, 1990.

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

Product Revision Date: 1/2/2009