

Monoclonal Anti-Poly Histidine Antibody

Product Number 00-1711 Clone Number AD1.1.10 Lot Number XXXX 100 ug Size Concentration 1 mg/ml

Formulation 0.2 µm filtered solution in 100 mM NaPi, 100 mM NaCl, 2 mM NaN₃

Storage

Specificity polypeptide containing a polyhistidine tag

Immunogen polyhistidine tagged polypeptide

IgG class mouse IgG₁

Preparation

This antibody was produced from a mouse hybridoma elicited from a mouse immunized with a polypeptide containing polyhistidine tag. The IgG fraction of ascites fluid was purified by Protein G affinity chromatography. The antibody pool was then purified with size exclusion chromatography (Superdex-200) to remove aggregated IgG.

Form/Shipping & Storage

Supplied as a 0.2 µm filtered solution in 100 mM sodium phosphate, 100 mM sodium chloride and 2 mM sodium azide. Can be stored at 2-8°C for at least three months without detectable loss of activity.

<u>Specificity</u>
This antibody has been selected for its ability to detect proteins containing accessible consecutive histidine regions. The antibody has been shown to detect polyhistidine tags localized at both the amino- or carboxyl-terminus.

For technical inquiries: info@columbiabiosciences.com For sales inquiries: sales@columbiabiosciences.com