



## Monoclonal Anti-Poly Histidine Antibody

Product Number	FITC-1711
Clone Number	AD1.1.10
Conjugate	FITC
Lot Number	XXXX
Size	100 ug
Concentration	0.1 mg/mL
Formulation	0.2 µm filtered solution in PBS, pH7.4 with 1% BSA
Storage	4-8°C or -20°C, store in the dark.
Specificity	polypeptide containing a polyhistidine tag
Immunogen	polyhistidine tagged polypeptide (PAX6 transcription factor linked to histidine tag)
IgG class	mouse IgG <sub>1</sub>
Applications	Flow Cytometry (1/10 dilution), Immunofluorescence (assay dependent) Western Blot (1:250 dilution).

### Preparation

This antibody was produced from cell culture elicited from a mouse immunized with a polypeptide containing polyhistidine tag. The cell culture was purified on Protein G and the resulting antibody pool was then polished by size exclusion chromatography (Superdex-200) to remove aggregated IgG.

### Form/Shipping & Storage

Supplied as a 0.2 µm filtered solution in PBS, pH 7.4 with 1%BSA. Can be stored at 2-8°C.

### Specificity

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.

4985 Winchester Blvd. Frederick, MD 21703  
301.732.5415

[www.columbiabiosciences.com](http://www.columbiabiosciences.com)

For technical inquiries: [info@columbiabiosciences.com](mailto:info@columbiabiosciences.com)

For sales inquiries: [sales@columbiabiosciences.com](mailto:sales@columbiabiosciences.com)