
Profinia™ His Antibody

Catalog Number
620-0203

BIO-RAD

Specifications

Source	Affinity-purified mouse IgG2b
Formulation	Supplied at 1 mg/ml in phosphate buffered saline, pH 7.4 with 0.05% NaN ₃ .
Specificity	The His-tag monoclonal antibody recognizes overexpressed 6x His fusion proteins.
Storage and stability	The antibody is shipped on blue ice and should be stored at -20°C for long-term storage. The antibody is stable for at least 1 year when stored at -20°C (see label for exact expiration). To avoid freeze-thaw cycles, aliquot the antibody. Working solutions can be stored at 4°C for up to 3 months.

Application

For colorimetric western blotting applications, use at 1:1,000–1:5,000 dilution. For chemiluminescent western blotting applications, use at 1:5,000 dilution.

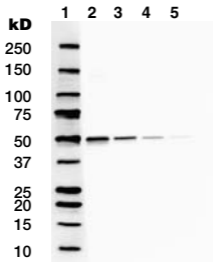


Fig. 1. Chemiluminescent detection of a 51 kD His tagged protein using the anti-His antibody at a 1:5,000 dilution. Lane 1: Precision Plus Protein™ standards detected with StrepTactin conjugate, Lanes 2–5: 250 ng, 60 ng, 15 ng, and 4 ng of the 51 kD target protein.

StrepTactin is a trademark of Institut für Bioanalytik GmbH.