
**Profinity eXact™ Tag
Antibody**

**Catalog Number
156-3004**

BIO-RAD

Specifications

Source	Affinity-purified mouse IgG1.
Formulation	Supplied at 1 mg/ml in phosphate buffered saline, pH 7.4 with 0.05% ProClin950.
Specificity	The Profinity eXact™-tag monoclonal antibody recognizes the 8.2 kD eXact tag domain in Profinity eXact-tagged 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). Aliquot the antibody to avoid freeze-thaw cycles. Working solutions can be stored at 4°C for up to 3 months.

Application

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

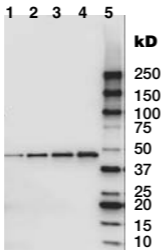


Fig. 1. Chemiluminescent detection of a 40 kD eXact-MBP fusion protein using the anti-eXact monoclonal antibody at 1:2,000 dilution.

Lanes 1–4: 30 ng, 60 ng, 100 ng, and 150 ng of the 40 kD fusion protein; Lane 5: Precision Plus Protein™ WesternC standards detected with StrepTactin conjugate.

StrepTactin is a trademark of Institut für Bioanalytik GmbH.

