

Results

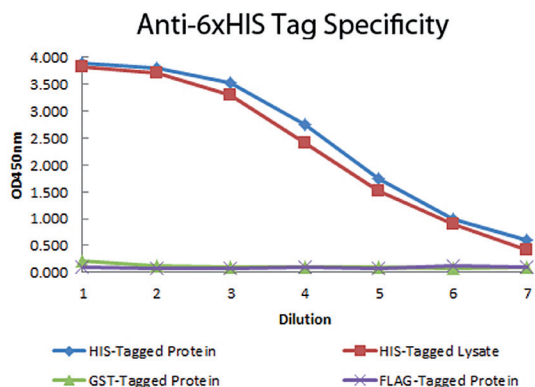


Figure 1. 6X His Tag antibody ELISA results

ELISA using the monoclonal 6xHis Tag Antibody. Antigen: HIS-tagged purified protein and E. coli cell lysates expressing HIS-Tagged construct, GST- and FLAG-tagged purified proteins. Coating amount: 0.15ug per well. Primary antibody: 6xHis Tag antibody at 100ug/mL. Dilution series: 2-fold. Mid-point concentration: 200ng/mL. Secondary antibody: Peroxidase mouse secondary antibody at 1:10,000.

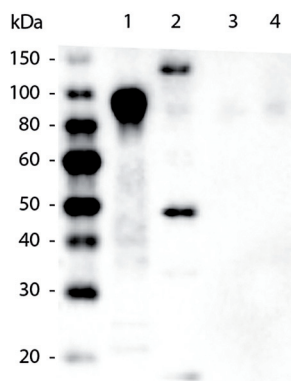


Figure 2. 6X His Tag antibody western blot results

Western Blot of Mouse anti-6xHis Tag Antibody. Lane 1: 100ng Purified histidine-tagged recombinant protein. Lane 2: 200ng E. coli cell lysate containing histidine-tagged expression construct. Lane 3: 100ng Purified GST-tagged recombinant protein. Lane 4: 100ng Purified FLAG-tagged recombinant protein. Primary antibody used at a 1:5,000 dilution overnight at 4°C. Secondary antibody: Peroxidase mouse secondary antibody at 1:20,000 for 30 min at RT.

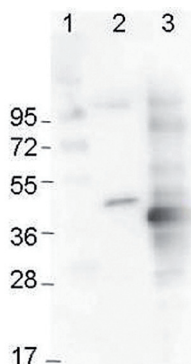


Figure 3. 6X His Tag antibody western blot results

Western Blot using the monoclonal 6x-His Epitope Tag Antibody showing detection of the 6xHis sequence on N-terminally-tagged (lane 2) and C-terminally-tagged recombinant proteins (lane 3). A molecular weight marker is shown in lane 1.

Diagenode sa. BELGIUM | EUROPE

LIEGE SCIENCE PARK
Rue Bois Saint-Jean, 3
4102 Seraing (Ougrée) - Belgium
Tel: +32 4 364 20 50
Fax: +32 4 364 20 51
orders@diagenode.com
info@diagenode.com

Diagenode Inc. USA | NORTH AMERICA

400 Morris Avenue, Suite 101
Denville, NJ 07834 - USA
Tel: +1 862 209-4680
Fax: +1 862 209-4681
orders.na@diagenode.com
info.na@diagenode.com

Last update: September 10, 2014