

Blue ladder monoclonal antibody

Cat. No. C15200213

Type: Monoclonal

Source: Mouse

Isotype: IgG1

Lot #: 003D

Size: 50 µl

Concentration: 2 µg/µl

Specificity: The antibody is specific for the blue colour used in protein MW markers. It does not react with human, mouse, rat, chicken, hamster, monkey, yeast and E. coli proteins.

Purity: Protein A purified monoclonal antibody in PBS containing 0.05% azide.

Storage: Store at -20°C; for long storage, store at -80°C. Avoid multiple freeze-thaw cycles.

Precautions: This product is for research use only. Not for use in diagnostic or therapeutic procedures.

Description: Monoclonal antibody raised in mouse against the blue colour used in many prestained protein MW markers. This antibody can be used to visualize the marker bands after WB.

Applications

	Suggested dilution/amount	References
Western blotting	1:1,000 - 1:5,000	Fig 1

Target description

Most prestained protein MW markers used in PAGE contain fragments that are labelled with a blue dye. This antibody specifically reacts with this blue dye, enabling to directly visualize the different marker fragments on the blot. This makes the positioning of the marker on the blot significantly more easy and accurate.

Results

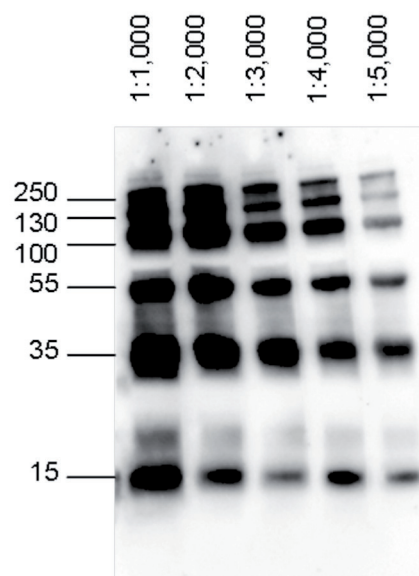


Figure 1. Western blot analysis using the Diagenode Blue ladder monoclonal antibody

5 μ l of the Page Ruler Plus Prestained Protein Ladder (Thermo Scientific, cat. No. 26619) were loaded on a gel and analyzed by Western blot with the Diagenode Blue ladder antibody (Cat. No. C1520213) used at different dilutions. Figure 1 shows that even with a 1:5,000 dilution all blue prestained bands of the molecular weight marker are clearly visible after a 30" exposure, a typical exposure time for WB analysis.

Western Blot protocol

1. Perform SDS-PAGE with a blue pre-stained protein molecular weight marker of your vendor of choice and your samples. Transfer proteins to a PVDF membrane.
2. Block the membrane in TBS-T (TBS + 0.05% Tween-20) containing 5% milk powder for 1h at room temperature.
3. Incubate the membrane overnight at 4°C with your primary antibody diluted in TBS-T + 5% milk. Incubate the lane with the molecular weight marker with the Blue ladder antibody diluted 1:1,000 to 1:5,000 in TBS-T + 5% milk. Alternatively, the complete membrane can be incubated with both antibodies together.
4. Wash the membrane and the lane with the molecular weight marker 3 times 10 minutes with TBS-T + 5% milk at room temperature.
5. Incubate the membrane and the lane with the molecular weight with a secondary HRP-coupled antibody diluted 1:10,000 to 1:20,000 in TBS-T + 5% milk for 1 hour at room temperature.
6. Wash the membrane and the lane with the molecular weight marker 3 times 10 minutes with TBS-T at room temperature.
7. Visualize your protein together with the blue pre-stained protein marker bands by enhanced chemiluminescence (ECL).

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: January 13, 2015