

PRODUCT DESCRIPTION

VisUBlock™ Mouse on Mouse Blocking Reagent effectively blocks endogenous IgG, which enables staining of mouse tissue with mouse IgG primary antibodies and eliminates unwanted background staining. Endogenous IgG poses a significant issue when using mouse IgG primary antibodies on mouse tissue: anti-mouse secondary detection antibodies will not only bind to mouse IgG primary antibodies, but also to endogenous tissue IgG. Therefore, if endogenous IgG is not blocked prior to incubation with secondary detection antibodies, intense background staining occurs which can mask specific staining generated by the primary antibody. 1.0 mL of concentrate is enough to stain 100-250 tissue sections when diluted.

INTENDED USE

VisUBlock™ Mouse on Mouse Blocking Reagent is for use with primary antibodies from a mouse host and mouse tissue.

STABILITY & STORAGE

Upon receipt, store at 2-8 °C for up to 12 months.

TECHNICAL HINTS & LIMITATIONS

- FOR LABORATORY RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES.
- Results may vary due to variations in tissue staining conditions.

REAGENT PREPARATION

Mix 30 µL of VisUBlock™ Mouse on Mouse Blocking Reagent with 1.0 mL of preferred buffer such as PBS or TBS. This will produce approximately 1.0 mL of working solution.

PROTOCOL

1. Prior to primary antibody incubation, apply Mouse on Mouse IgG Blocking Reagent to tissue section. Incubate 1 hour at room temperature.
2. Tap off, do not rinse, and continue with primary antibody incubation and detection antibody incubation according to protocol set in lab.