

BLOXALL® Endogenous Blocking Solution

Peroxidase and Alkaline Phosphatase

Cat. No. SP-6000

Storage Refrigerate 2-8 °C

Description BLOXALL Endogenous Peroxidase and Alkaline Phosphatase Blocking Solution is a dual enzyme blocking reagent that inactivates endogenous peroxidase, pseudoperoxidase, and alkaline phosphatase in formalin-fixed, paraffin-embedded tissue sections, frozen tissue sections, and cell preparations. This reagent is provided as a ready-to-use reagent in a convenient dropper bottle.

Components

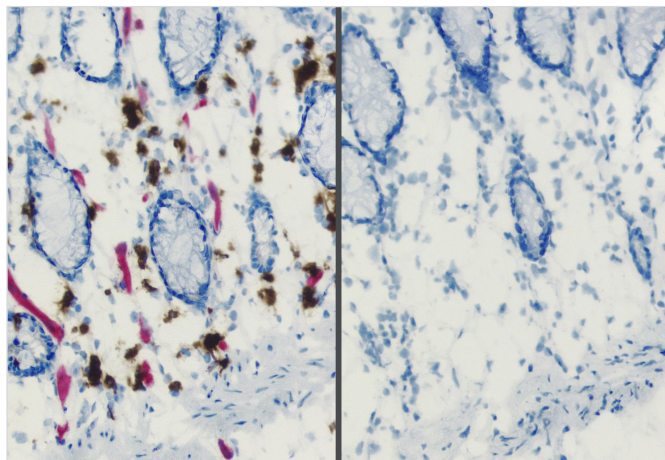
Product Name	Volume
BLOXALL Endogenous Blocking Solution	100 ml

Guidelines for Use

Prior to the application of primary antibody, apply BLOXALL Endogenous Blocking Solution to tissue sections or cell preparations and incubate for 10 minutes at room temperature. Wash sections in an appropriate wash buffer such as Tris buffer, PBS, or TBS for 5 minutes. Incubation time can be shortened based on individual application.

Appropriate care should be exercised when using this reagent to avoid contact with eyes and skin. Dispose in accordance with local regulations.

Detailed product listings, specifications, protocols and additional information is available on our website: vectorlabs.com



Endogenous alkaline phosphatase (AP) and peroxidase (HRP) activities in frozen intestine revealed with Vector® Red AP Substrate, red and ImmPACT® DAB HRP Substrate, brown (left). Same substrates used on BLOXALL Solution-treated issue (right).