



CERTIFICATE OF ANALYSIS

Product **BLOXALL™ Endogenous Peroxidase and Alkaline Phosphatase Blocking Solution**

Catalog No. SP-6000

Amount 100 ml

Lot No. X1010

Storage Refrigerate 4° C - 8° C

BLOXALL™ Blocking Solution will inactivate endogenous peroxidase, pseudoperoxidase, and alkaline phosphatase in formalin-fixed, paraffin-embedded tissue sections, frozen tissue sections, and cell preparations.

BLOXALL™ Blocking Solution is provided ready-to-use in a dropper bottle.

Prior to the application of primary antibody, apply BLOXALL™ 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.