

VitroView™ Antibody Dilution Buffer (100 ml)

Cat. No. VB-6002

Introduction

This antibody dilution buffer is ready-to-use solution for dilution of primary and secondary antibodies in immunoassays including immunofluorescence (IF), Immunohistochemistry (IHC) and Western blot (WB).

Kit Components

BSA (IHC grade, IgG-Free, Protease -free)
cold fish skin gelatin
Sodium azide
0.01M PBS, pH7.2

Storage

Store at 4°C.

Application

IHC, ICC/IF, WB. The usefulness of this product in other applications has not been determined.

Recommended Assay Procedure for Immunohistochemistry

1. After DAB staining, incubate sections with 3 drops of RTU hematoxylin solution for 1-2 minutes
2. Rinse in tap water 2x2 min.
3. Dehydrate through 75% ethanol for 2 min, 95% ethanol for 2 min, and 100% ethanol for 2x3min. Clear in xylene for 2x5min.
4. Coverslip with mounting medium.

Note

1. The sodium azide is an inhibitor of HRP. Do not use it to dilute HRP conjugated antibody.
2. The phosphate in PBS is an inhibitor of Alkaline Phosphatase (AP). Do not use it to dilute AP conjugated antibody.

Precautions

Handle with care. Avoid contact with eyes, skin and clothing. Do not ingest. Wear gloves.