VectaPlex™ Antibody Removal Kit (VRK-1000)



Together we breakthrough™

Can I use VectaPlex™ on frozen sections?

No. During development, VectaPlex $^{\text{TM}}$ was optimized for FFPE tissues, where the power to remove antibodies at room temperature was balanced with the need to preserve tissue morphology. For frozen tissues, VectaPlex $^{\text{TM}}$ will remove the tissue from the slide.

Can VectaPlex™ remove antibodies in immunohistochemistry (IHC) applications with chromogenic staining?

VectaPlex[™] Antibody Removal Kit is designed to be used in fluorescent workflows such as immunofluorescence or TSA. It has not been validated or confirmed for immunohistochemistry (IHC) applications. VectaPlex[™] may remove substrates already deposited on tissue samples.

I am using TrueView to quench autofluorescence, will TrueVIEW be stripped by VectaPlex™ reagents?

We used TrueView in our VectaPlexTM IF workflow to suppress autofluorescence. Application of VectaPlexTM removes some, but not all, of the TrueView reagent. Since some TrueView remained after VectaPlexTM application, we did not add TrueView in every cycle, but only as needed to suppress autofluorescence.

I accidentally put my VectaPlex™ kit in the freezer. Can I still use it?

Yes. VectaPlex[™]can withstand a freeze-thaw cycle and still perform as expected, but the recommended storage is at ambient temperature (15-25 °C). Additionally, the kit will still perform properly if placed at 4 °C.

Is there a maximum number of times you can use VectaPlex™ on one section?

VectaPlex[™] has been tested for up to 6 cycles with no ill effects. We don't expect there to be issues with increasing the number of cycles, but we suggest testing with controls for adequate comparisons. We recommend running experiments to confirm that the use of more cycles does not impact the antigenicity or tissue morphology.

Can VectaPlex™ be used on an autostainer?

We have not tested this product on automated systems and cannot guarantee instrument compatibility at this time.

Does VectaPlex™ work for all species of antibodies?

VectaPlex[™] has been tested with mouse and rabbit primaries, and with horse and goat secondaries. We expect it to work well with other species of antibodies as well.

Does VectaPlex™ work on different tissue types?

VectaPlex[™] has been tested on and works well with various FFPE tissues, including FFPE breast and lung tissue which are very fragile.

Is VectaPlex™ compatible with all antibodies?

VectaPlex[™] has been tested on more than 30 common antibodies and has worked on them without issue.

In an IF protocol, will the VectaPlex[™] Antibody Removal Kit also strip the autofluorescence from the tissue?

VectaPlex[™] does not appear to reduce autofluorescence background.

Is there a protocol for removal of coverslips between VectaPlex™ cycles?

We recommend using VectaShield® Plus Antifade Mounting Media with DAPI (H-2000), which is an aqueous non-hardening mounting medium. After each imaging cycle, the coverslip can be easily removed by incubating in PBS and the residual VectaShield® removed with a PBS wash.