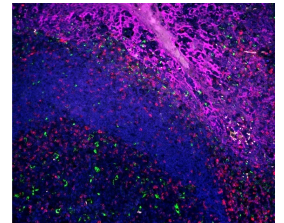
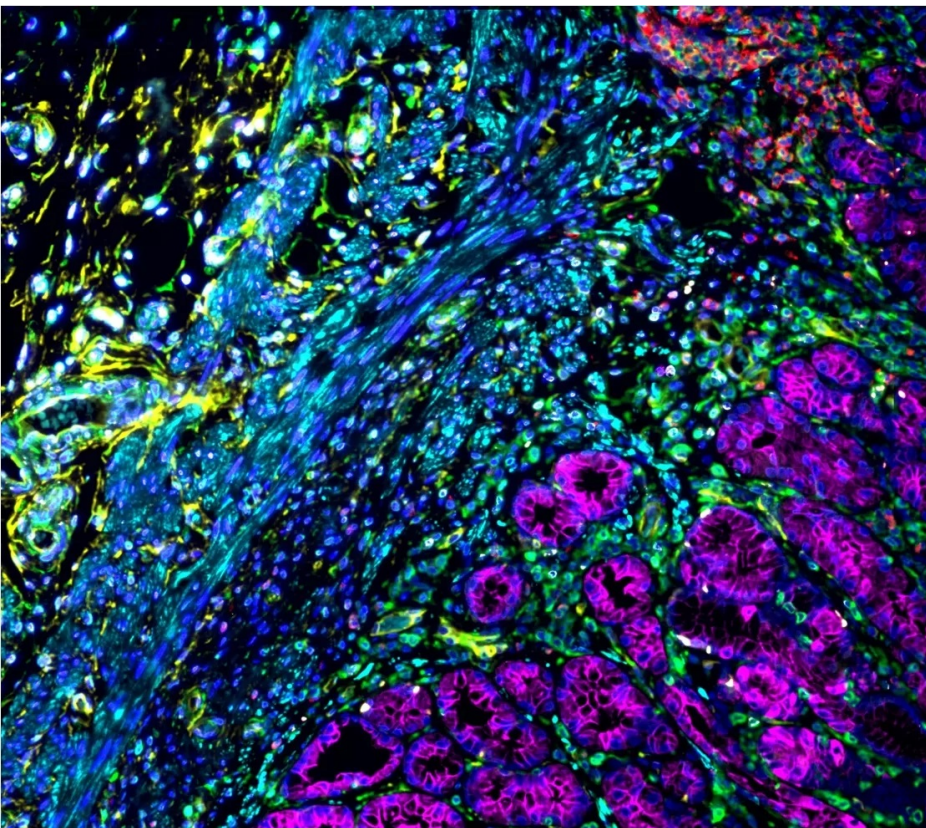




VectaPlex™ Antibody Removal Kit

VRK-1000

Product Images



Short Description

VectaPlex Antibody Removal kit is a ready-to-use kit designed to remove antibodies and other non-covalently bound detection reagents from formalin-fixed paraffin-embedded (FFPE) tissue sections in immunofluorescence or tyramide signal amplification staining workflows.

Additional Information

Unit Size	1 kit
Applications	Immunofluorescence, Tyramide Signal Amplification
Recommended Usage	(15–25° C)
Target Species	Universal
Format	Ready-to-Use

