

VECTOR LABORATORIES

Enzyme Substrate Combinations

This table is designed as a reference for optimal multiple labeling, since the order of labeling and the relative intensities of the two color precipitates may significantly affect the quality and labeling pattern of each antigen in the stained section.

Second Substrate \ First Substrate	Enzyme Substrate Combinations											
	Vector® Red (red)	Vector® Blue (blue)	Vector® Black (brown/black)	BCIP/NBT (blue/violet)	Vector® VIP (purple)	ImmPACT™ DAB, DAB (brown)	DAB-Ni (gray/black)	Vector® NovaRED™ (red)	Vector® SG (blue/gray)	AEC (red)	TMB (blue)	
Vector® Red (red) Cat. No. SK-5100		-	-	-	-	+	+	-	+	-	-	
Vector® Blue (blue) Cat. No. SK-5300	+		-	-	+	+	+	+	+	+	-	
Vector® Black (brown/black) Cat. No. Sk-5200	+	+		+	+	-	-	-	-	+	-	
BCIP/NBT (blue/violet) Cat. No. SK-5400	+	-	-		+	+	+	+	+	+	-	
Vector® VIP (purple) Cat. No. SK-4600	-	+	-	-		+	+	-	+	-	-	
ImmPACT™ DAB, DAB (brown) Cat. Nos. SK-4105, SK-4100	+	+	-	+	+		-	-	+	+	+	
DAB-Ni (gray/black) Cat. No. SK-4100	+	-	-	-	+	+		+	-	+	+	
Vector® NovaRED™ (red) Cat. No. SK-4800	-	+	-	+	-	+	+		+	-	+	
Vector® SG (blue/gray) Cat. No. Sk-4700	+	-	-	-	+	+	-	-		+	-	
AEC (red) Cat. No. SK-4200	-	-	-	-	-	+	-	-	+		-	
TMB* (blue) Cat. No. SK-4400	-	-	-	-	-	-	-	-	-	-		

KEY:

Alkaline Phosphatase Peroxidase

+ indicates good contrast, - indicates incompatibility of substrates for various reasons

* TMB does not provide optimal reactivity in double label applications.