

VECTOR LABORATORIES

Enzyme Substrate Combinations

This table shows the substrates that are recommended for multiple antigen labeling and the order in which they are optimally used.

Second Substrate \ First Substrate	Vector® Red (magenta) SK-5100	Vector® Blue (blue) SK-5300	BCIP/NBT (indigo) SK-5400	VIP/ ImmPACT™ VIP (purple) SK-4605, SK-4600	DAB/ ImmPACT™ DAB (brown) SK-4105, SK-4100	DAB-Ni (gray-black) SK-4100	NovaRED™/ ImmPACT™ NovaRED™ (red) SK-4805, SK-4800	SG/ ImmPACT™ SG (blue-gray) SK-4705, SK-4700	AEC/ ImmPACT™ AEC (red) SK-4205, SK-4200	ImmPACT™ AMEC <i>Red</i> (red) SK-4285
Vector® Red (magenta) Cat. No. SK-5100		–	–	–	+	+	–	+	–	–
Vector® Blue (blue) Cat. No. SK-5300	+		–	+	+	+	+	+	+	+
BCIP/NBT (indigo) Cat. No. SK-5400	+	–		+	+	+	+	+	+	+
VIP/ImmPACT™ VIP (purple) Cat. No. SK-4605, SK-4600	–	+	–		+	+	–	+	–	–
DAB/ImmPACT™ DAB (brown) Cat. No. SK-4105, SK-4100	+	+	+	+		–	–	+	+	+
DAB-Ni (gray-black) Cat. No. SK-4100	+	–	–	+	+		+	–	–	–
NovaRED™/ImmPACT™ NovaRED™ (red) Cat. No. SK-4805, SK-4800	–	+	+	–	+	+		+	–	–
SG/ImmPACT™ SG (blue-gray) Cat. No. SK-4705, SK-4700	+	–	–	+	+	–	–		+	+
AEC/ImmPACT™ AEC (red) Cat. No. SK-4205, SK-4200	–	–	–	–	+	–	–	+		–
ImmPACT™ AMEC <i>Red</i> (red) SK-4285	–	–	–	–	+	–	–	+	–	

Key:

Alkaline Phosphatase

Peroxidase

+ Indicates good contrast

– Indicates incompatibility of substrates for various reasons