

间接 ELISA

abcam[®]
discover more

Coat plate with antigen
2 hr RT or 4°C overnight



Wash plates with PBS 0.2% Tween 20 4 times

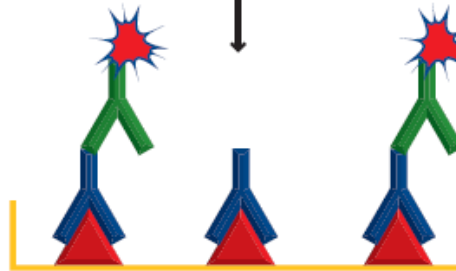
Block with 5% serum or BSA 2 hr or overnight 4°C

Incubate with
primary antibody
2 hr RT or 4°C overnight



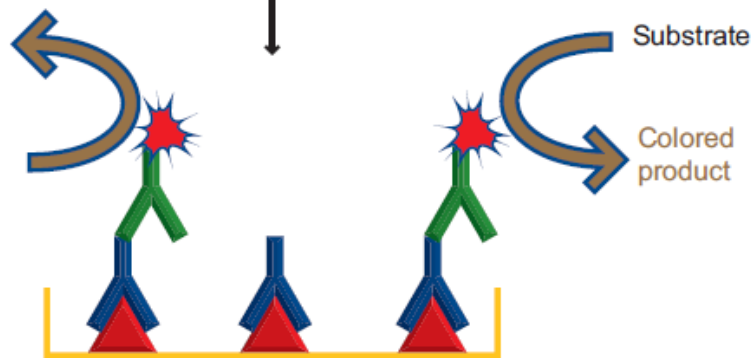
Wash plates with PBS 0.2% Tween 20 4 times

Add conjugated
secondary antibody
Incubate 1 - 2 hr



Wash plates with PBS 0.2% Tween 20 4 times

Enzymatic detection
Follow manufacturers
recommendations



Read absorbance on ELISA plate
reader and analyze results

Key

