

Sandwich ELISA

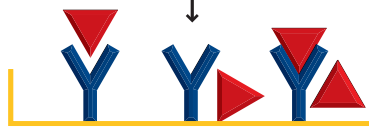
Incubate with coating antibody in bicarbonate buffer overnight 4°C



Wash plates with PBS 0.2% Tween 20 2 times

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

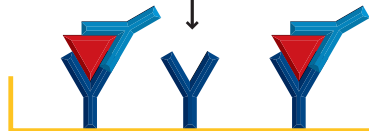
Add sample to wells and dilute down the plate



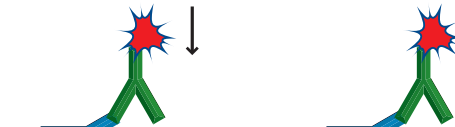
Wash plates with PBS 0.2% Tween 20 2 times



Incubate with detection antibody 2 hr RT



Wash with PBS 0.2% Tween 20 4 times



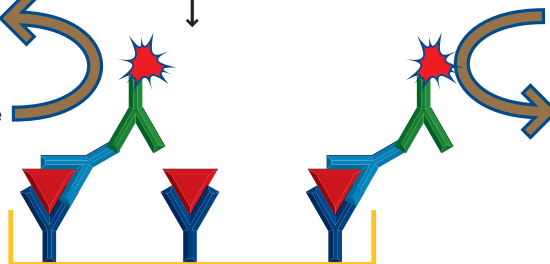
Incubate with conjugated secondary antibody 30 min to 2 hr RT



Wash plates with PBS 0.2% Tween 20 4 times

Colored product





Substrate



Enzymatic detection
Follow manufacturers recommendations

Read absorbance on ELISA plate reader and analyze results

Key

-  Antigen
-  Primary capture antibody
-  Primary detection antibody
-  Conjugated Secondary antibody