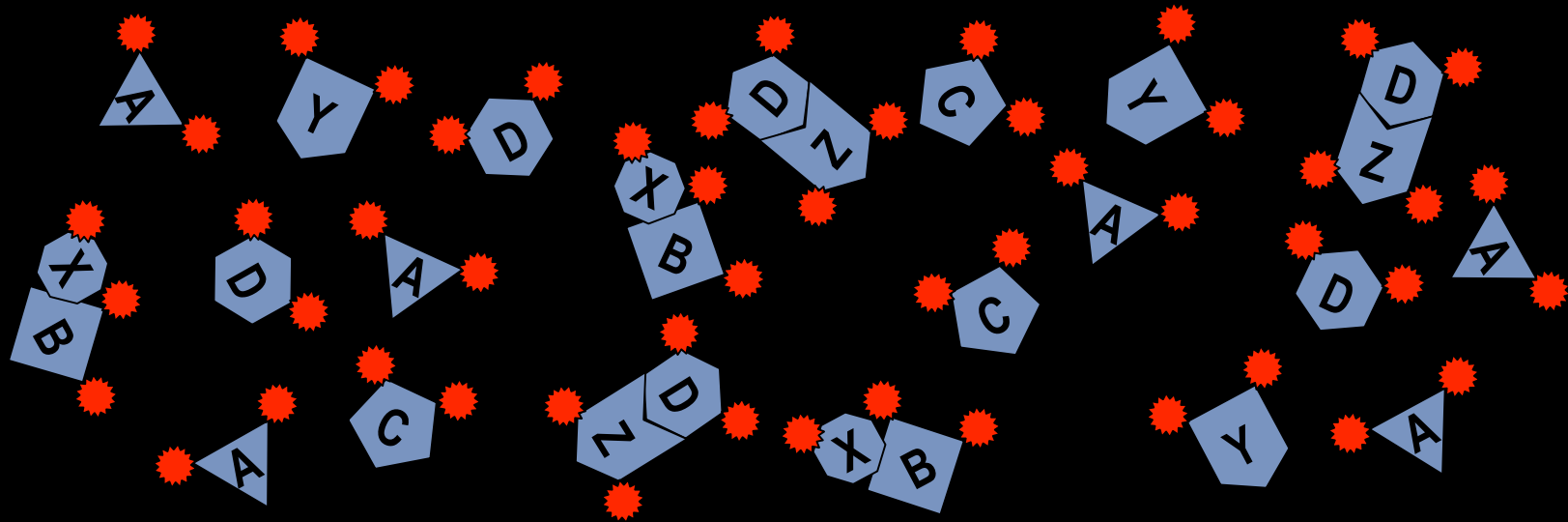


Kinex™ Antibody Microarrays

Step 2



- Label cell lysate proteins with fluorescent dye