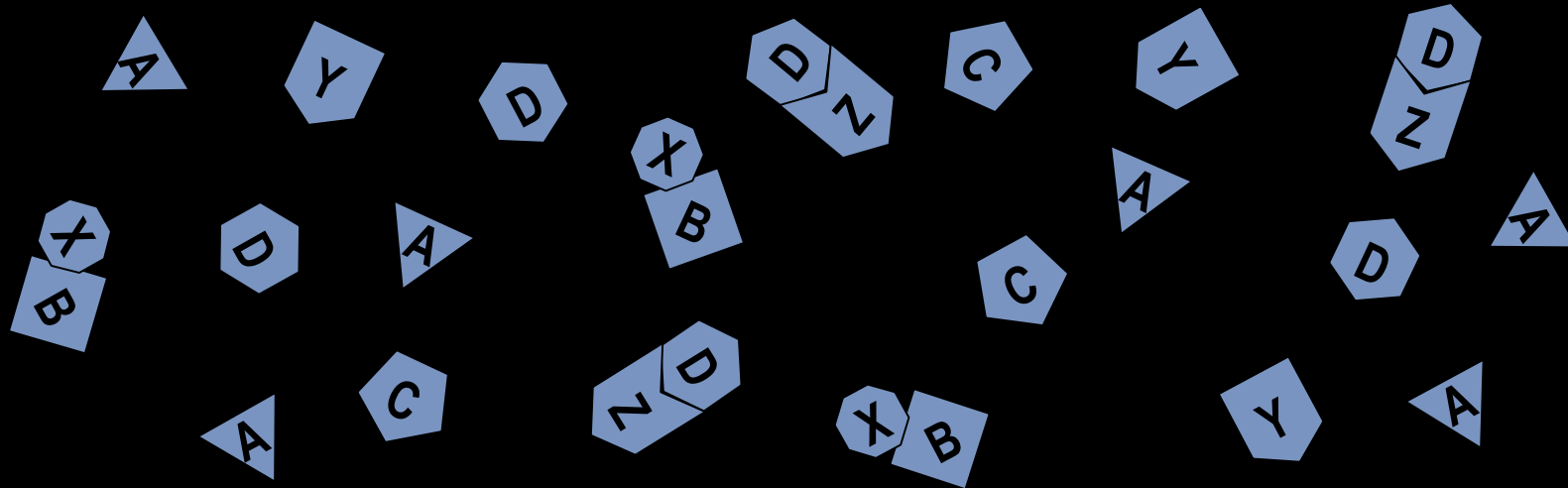


Kinex™ Antibody Microarrays

Step 1



- Start with cell lysate with thousands of different proteins

KINEXUS



www.kinexus.ca