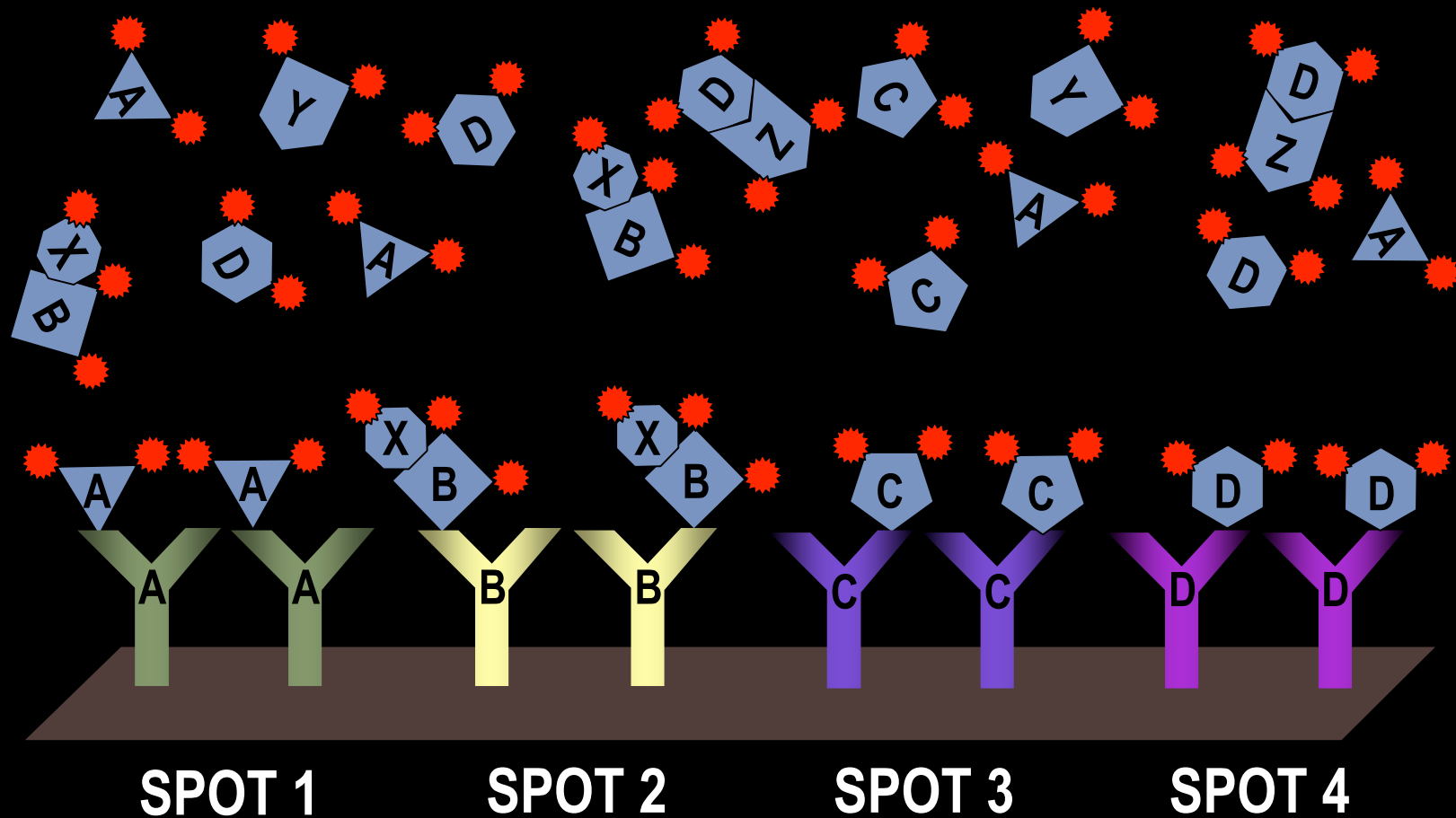


Kinex™ Antibody Microarrays

Step 4



KINEXUS



www.kinexus.ca