

Reverse Microarray Format

 **Biotin**

 **Anti-phospho-site Antibody**

 **Dye**

 **Streptavidin**

- Biotinylated antibody for a specific target protein or phospho-site is used to probe an array of tissue lysates

