

Pan-Phospho Antibody in Sandwich Microarray Format



■ Antibody captured proteins are probed with a pan-specific phospho-site reporter antibody

