

Kinetworks™ Phospho-Site Multi-Immunoblot

Jurkat T Cells -Staurosporin Treated Overnight

