

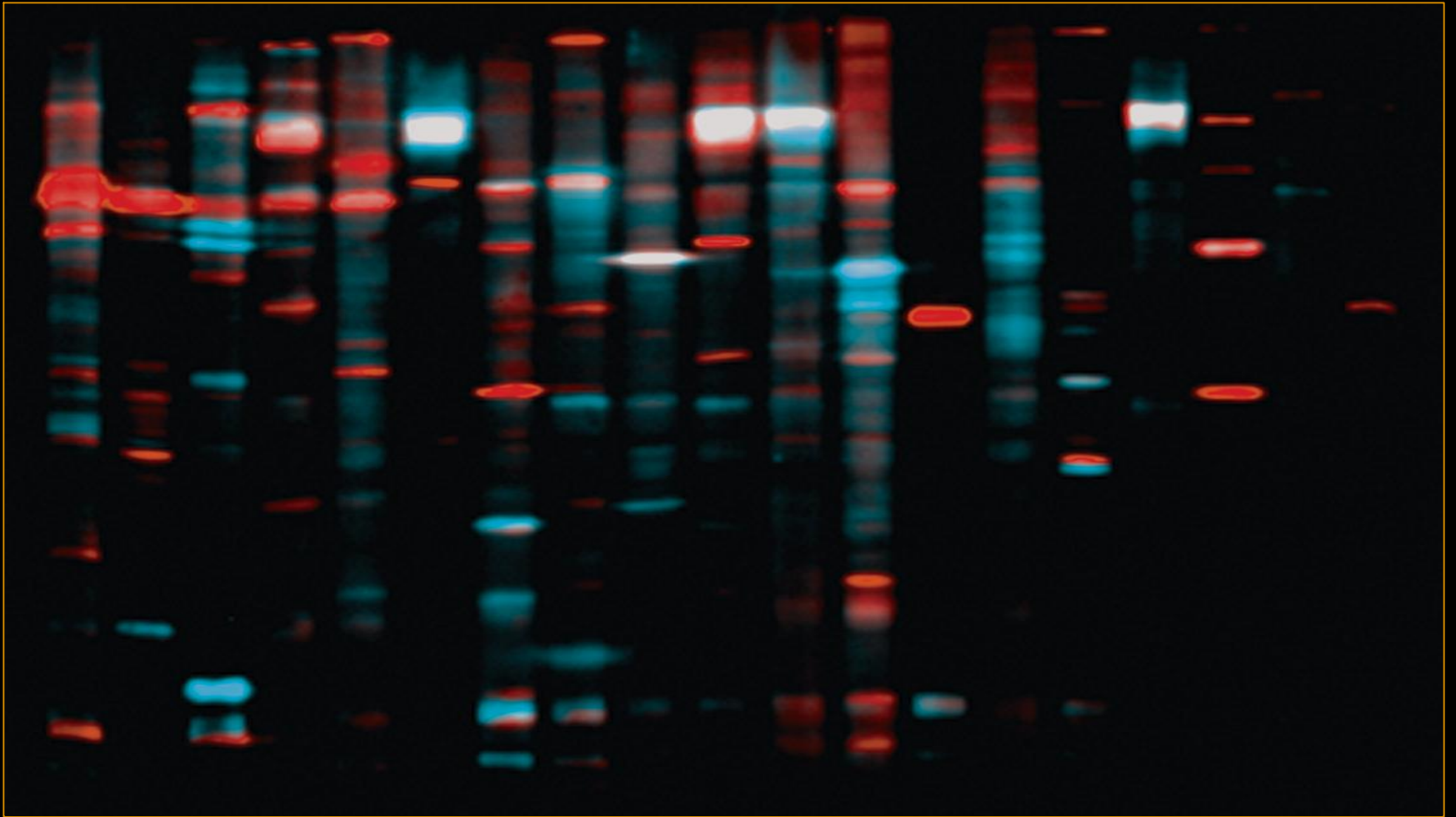
Human Jurkat T cells



Untreated 

KINEXUS

Kinetworks™ KPSS 8.0 Phospho-Site Screen



Human Jurkat T cells



Untreated  Staurosporin  Overlap 

KINEXUS

Kinetworks™ KPSS 8.0 Phospho-Site Screen



Human Jurkat T cells

Untreated ■ Staurosporin ■ Overlap ■

KINEXUS

Kinetworks™ KPSS 8.0 Phospho-Site Screen