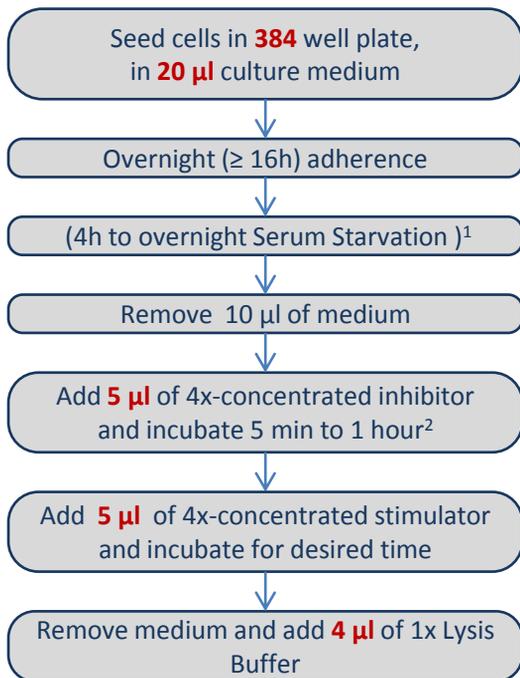
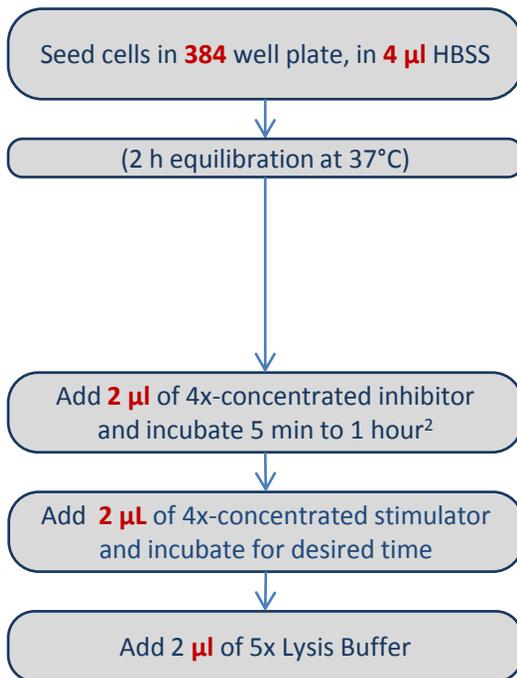


AlphaScreen® SureFire®: 1-plate/ 2-steps assay flowchart

Adherent Cells



Suspension Cells

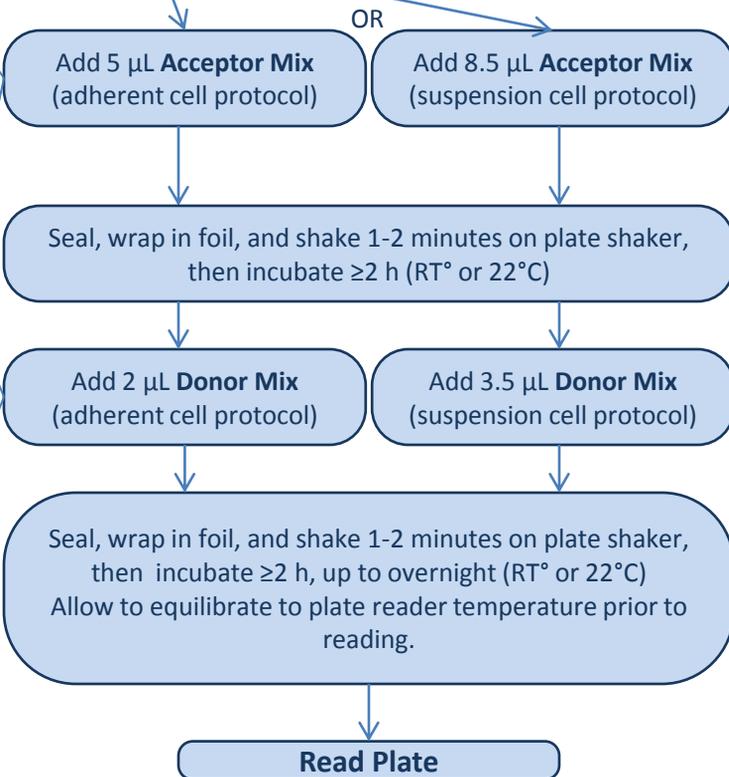


Incubate for 10 min on plate shaker (~350 rpm).³

In control wells, add positive or negative control lysates.

Acceptor Mix:	Typical volume ⁴	My volumes
Reaction Buffer	392 µL	
Activation Buffer	98 µL	
Acceptor Beads	10 µL	

Donor Mix:	Typical volume ⁴	My volumes
Dilution Buffer	190 µL	
Donor Beads	10 µL	



⁴ Important note: the volumes shown here are the ones for most common kits. For a few kits (e.g. c-Jun, cofilin, GSK3β, H3, MEK-1 and VEGFR2), the dilution factors are different. In this case please refer to the manual of the kits for volumes to be used.

¹ Depending on cell type and pathway analyzed.

² Depending on type of inhibitor used: 5 min is generally enough for receptor antagonists; more time is needed to block intracellular targets.

³ May stop and freeze lysates at -20°C if desired. If doing this, re-shake after thawing to ensure homogeneity of solution.