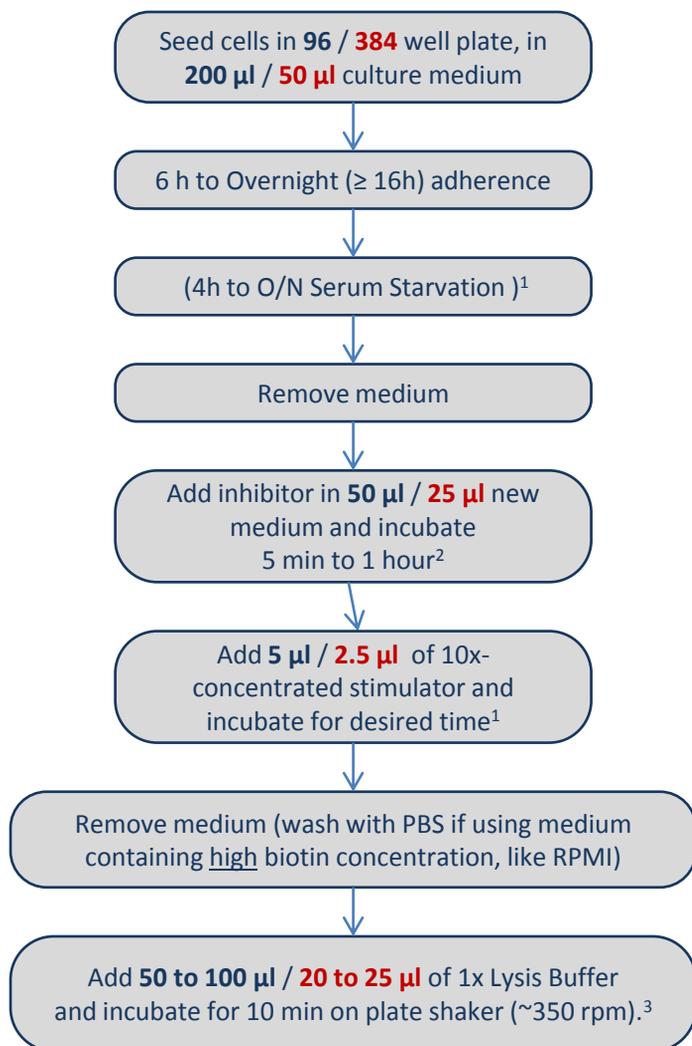
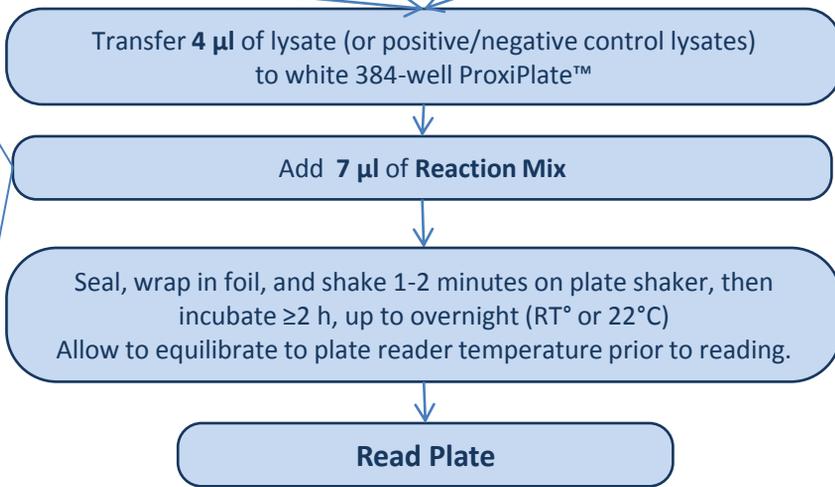
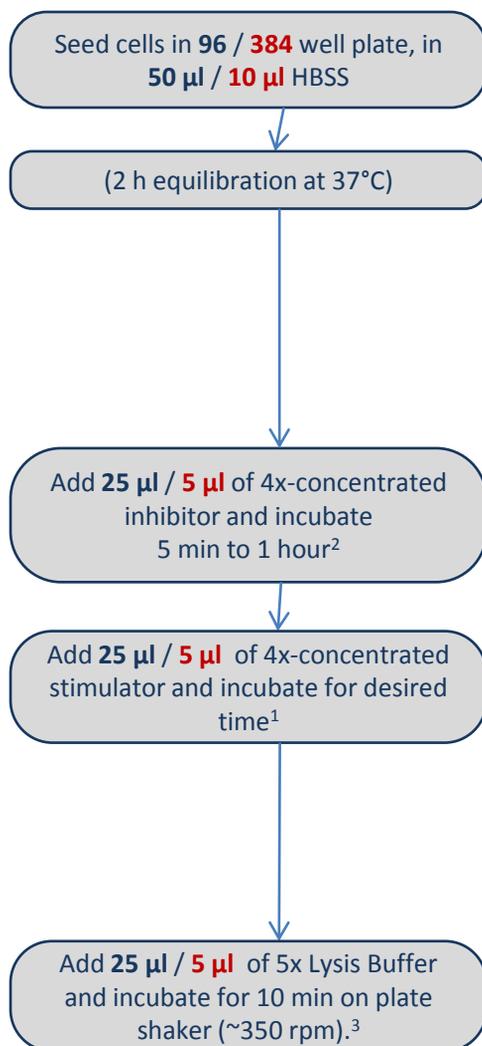


AlphaScreen® SureFire®: 2-plates/ 1-step assay flowchart (p-ERK1/2 and p-MEK1 kits only)

Adherent Cells



Suspension Cells



Reaction Mix:	Typical volume for p-ERK1/2 ⁴	Typical volume for p-MEK1 ⁴
Reaction Buffer	587 µL	636 µL
Activation Buffer	98 µL	49 µL
Acceptor Beads	10 µL	10 µL
Donor Beads	5 µL	5 µL

⁴ Important note: only the p-ERK and p-MEK1 assays can be run using a single step immunoassay. For other kits; a separate Acceptor Mix and Donor Mix procedure needs to be performed.

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.