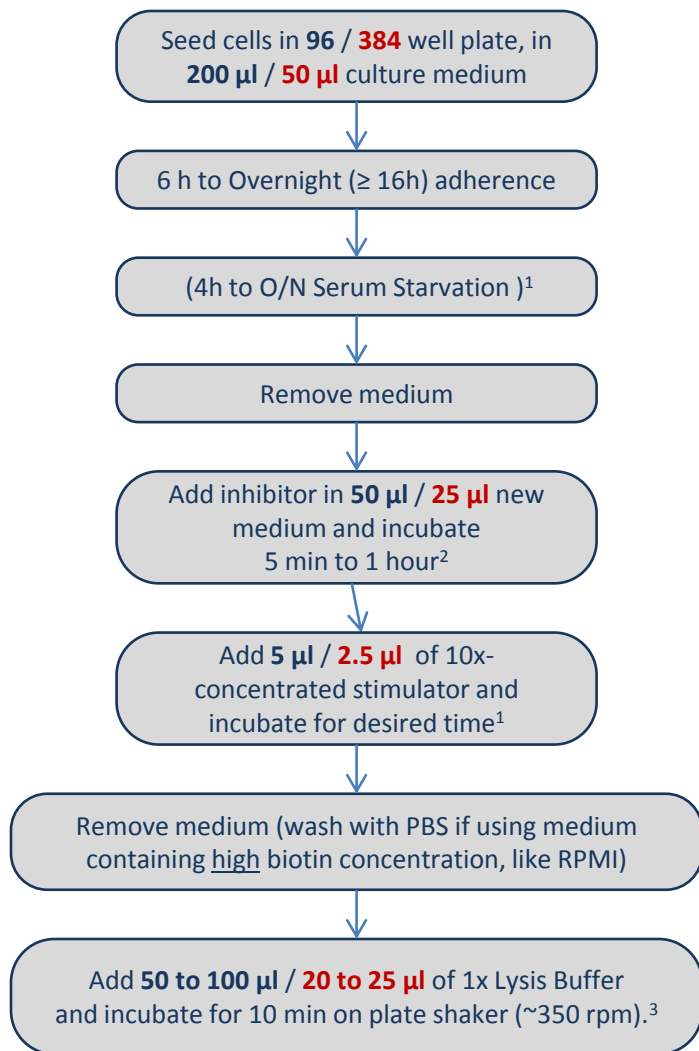
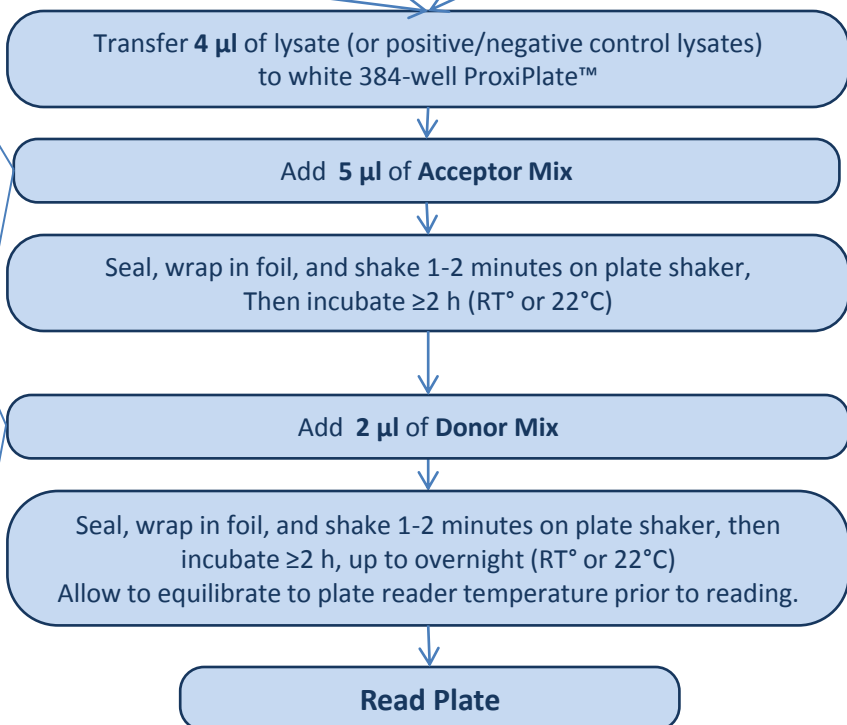
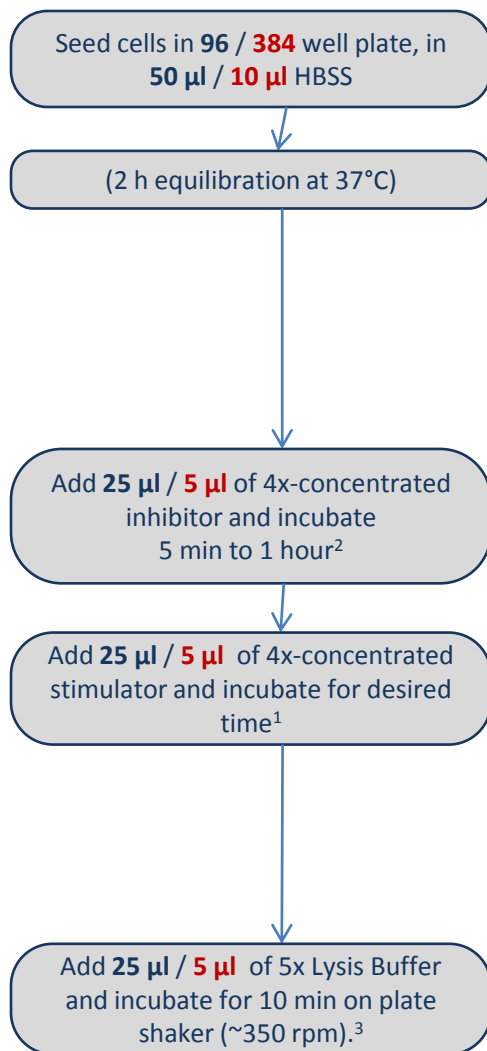


AlphaScreen® SureFire®: 2-plates/ 2-steps assay flowchart

Adherent Cells



Suspension Cells



| 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.