

## Flow Cytometry Protocol (clone 28-8, Ab205921)

### Flow Cytometry Materials

- Primary Antibody: Anti-PDL1 (Clone 28-8)
- Isotype Control: Rabbit IgG ([ab172730](#))
- Secondary Antibody: Donkey anti-Rabbit (Fab)'2-APC
- FACS Buffer: 1% BSA / 0.1% NaAzide in PBS (buffer)
- 96 well u-bottom plates, plate sealers, 50mL conical Tubes
- Human Gamma-Globulin
- L2987 (PD-L1 High) and A549 (PD-L1 Low) human tumor cells

### Cell Preparation

- L2987 and HT29 cells are harvested using 0.25% Trypsin
- Cells are counted, then centrifuge @ 300G for 5 min
- Wash cells in 10mL buffer
- Bring cells to  $10 \times 10^6$  cells/mL in buffer

### Staining

- Block Fc-receptors using 20 $\mu$ L / mL human gamma-globulin
- Incubate 20 min, on ice
- Transfer 100 $\mu$ L ( $1 \times 10^6$  cells) per well
- Add 1 $\mu$ g / well anti-PDL1 or isotype
- Incubate plate 30 min, on ice, in dark
- Add 100 $\mu$ L buffer, centrifuge 300G for 5 min, decant
- Add 100 $\mu$ L buffer, centrifuge 300G for 5 min, decant
- Dilute Secondary mAb 1:66 (45 $\mu$ L stock + 2.95mL buffer)
- Add 100 $\mu$ L diluted secondary mAb to each well
- Incubate plate 20 min, on ice, in dark
- Add 100 $\mu$ L buffer, centrifuge 300G for 5 min, decant
- Add 100 $\mu$ L buffer, centrifuge 300G for 5 min, decant
- Add 100 $\mu$ L 1% formaldehyde, incubate 20 min on ice in dark
- Add 100 $\mu$ L buffer, centrifuge 300G for 5 min, decant
- Add 200 $\mu$ L buffer per well, seal plate and store in fridge
- Read samples on BD FACS CANTO