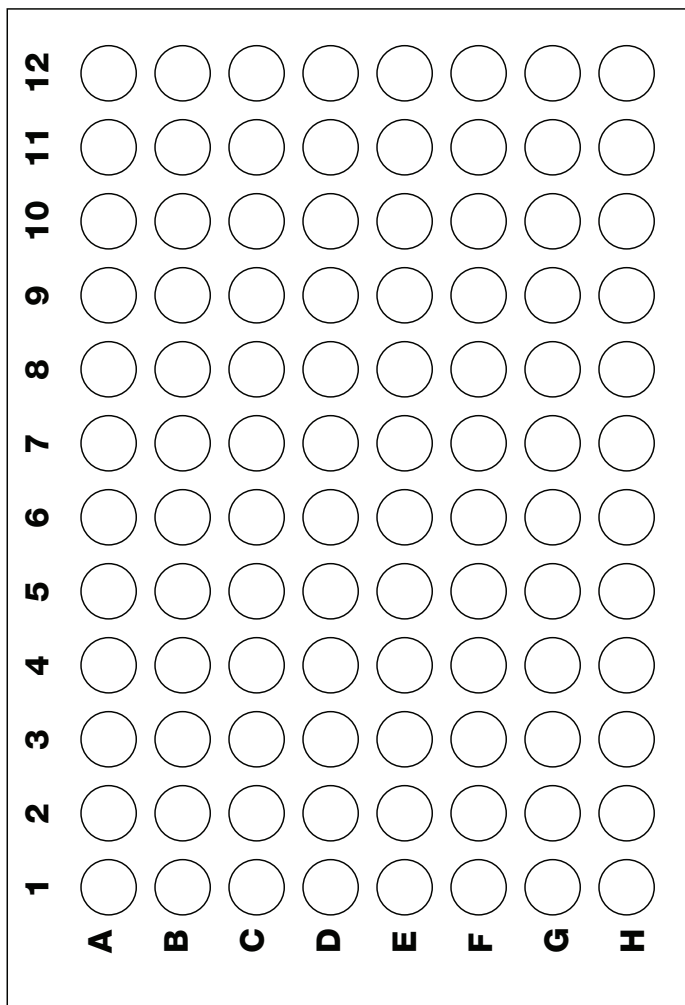


# Plate Layout Template



# Calculation Worksheet

## Plan Plate Layout

1. Fill out the 96-well plate template as instructed in the Plan Plate Layout section.

If using either the **Diabetes fixed panel (either human or mouse) or One Diabetes single set tube/analyte**, follow these directions:

Enter the number of wells that will be used in the assay: \_\_\_\_\_ (1)

## Calculations for Coupled Beads

1. Determine the volume of 1x coupled beads needed.

a) Each well requires 50  $\mu\text{l}$  of coupled beads (1x): \_\_\_\_\_ (1)  $\times$  50  $\mu\text{l}$  = \_\_\_\_\_  $\mu\text{l}$  (2)

b) Include a 20% excess to ensure enough volume: \_\_\_\_\_  $\mu\text{l}$  (2)  $\times$  0.20 = \_\_\_\_\_  $\mu\text{l}$  (3)

c) Total volume of 1x coupled beads: \_\_\_\_\_  $\mu\text{l}$  (2) + \_\_\_\_\_  $\mu\text{l}$  (3) = \_\_\_\_\_  $\mu\text{l}$  (4)

d) Volume of **20x coupled beads** stock: \_\_\_\_\_  $\mu\text{l}$  (4)/20 = \_\_\_\_\_  $\mu\text{l}$  (5)

e) Volume of **Assay Buffer** required: \_\_\_\_\_  $\mu\text{l}$  (4) \_\_\_\_\_  $\mu\text{l}$  (5) = \_\_\_\_\_ (6)

## Calculations for Detection Antibodies

2. Determine the volume of 1x detection antibody needed.

a) Each well requires 25  $\mu\text{l}$  of detection antibodies (1x): \_\_\_\_\_ (1)  $\times$  25  $\mu\text{l}$  = \_\_\_\_\_  $\mu\text{l}$  (7)

b) Include a 25% excess to ensure enough volume: \_\_\_\_\_  $\mu\text{l}$  (7)  $\times$  0.25 = \_\_\_\_\_  $\mu\text{l}$  (8)

c) Total volume of 1x detection antibodies: \_\_\_\_\_  $\mu\text{l}$  (7) + \_\_\_\_\_  $\mu\text{l}$  (8) = \_\_\_\_\_  $\mu\text{l}$  (9)

d) Volume of **20x Detection Antibodies** stock: \_\_\_\_\_  $\mu\text{l}$  (9)/20 = \_\_\_\_\_  $\mu\text{l}$  (10)

e) Volume of **Detection Antibody Diluent** required: \_\_\_\_\_  $\mu\text{l}$  (9) - \_\_\_\_\_  $\mu\text{l}$  (10) = \_\_\_\_\_  $\mu\text{l}$  (11)

## Calculations for Streptavidin-PE

3. Determine the volume of 1x streptavidin PE needed.

a) Each well requires 50  $\mu\text{l}$  of streptavidin PE (1x): \_\_\_\_\_ (1)  $\times$  50  $\mu\text{l}$  = \_\_\_\_\_  $\mu\text{l}$  (10)

b) Include a 25% excess to ensure enough volume: \_\_\_\_\_  $\mu\text{l}$  (10)  $\times$  0.25 = \_\_\_\_\_  $\mu\text{l}$  (11)

c) Total volume of 1x detection antibodies: \_\_\_\_\_  $\mu\text{l}$  (10) + \_\_\_\_\_  $\mu\text{l}$  (11) = \_\_\_\_\_  $\mu\text{l}$  (12)

d) Volume of **100x Streptavidin PE** required: \_\_\_\_\_  $\mu\text{l}$  (12) / 100 = \_\_\_\_\_  $\mu\text{l}$  (13)

e) Volume of **Assay Buffer** required: \_\_\_\_\_  $\mu\text{l}$  (12) \_\_\_\_\_  $\mu\text{l}$  (13) = \_\_\_\_\_  $\mu\text{l}$  (14)

If **multiplexing single set (singleplex) diabetes analytes**, follow these directions:

Enter the number of wells that will be used in the assay: \_\_\_\_\_ (1)

## Calculations for Coupled Beads

1. Determine the volume of 1x coupled beads needed.

a) Each well requires 50  $\mu\text{l}$  of coupled beads (1x): \_\_\_\_\_ (1)  $\times$  50  $\mu\text{l}$  = \_\_\_\_\_  $\mu\text{l}$  (2)

b) Include a 20% excess to ensure enough volume: \_\_\_\_\_  $\mu\text{l}$  (2)  $\times$  0.20 = \_\_\_\_\_  $\mu\text{l}$  (3)

c) Total volume of 1x coupled beads: \_\_\_\_\_  $\mu\text{l}$  (2) + \_\_\_\_\_  $\mu\text{l}$  (3) = \_\_\_\_\_  $\mu\text{l}$  (4)

d) Enter the number of diabetes single set (or analytes) tubes that will be multiplexed = \_\_\_\_\_(5)

e) Volume of **20x Coupled Beads** required from **each diabetes coupled beads tube**:

\_\_\_\_\_  $\mu\text{l}$  (4) /20 = \_\_\_\_\_ $\mu\text{l}$  (6)

f) Total volume of diabetes bead stock required: \_\_\_\_\_(5)  $\times$  \_\_\_\_\_  $\mu\text{l}$  (6) = \_\_\_\_\_  $\mu\text{l}$  (7)

g) Volume of **Assay Buffer** required: \_\_\_\_\_  $\mu\text{l}$  (4) \_\_\_\_\_  $\mu\text{l}$  (7) = \_\_\_\_\_  $\mu\text{l}$  (8)

## Calculations for Detection Antibodies

2. Determine the volume of 1x detection antibody needed.

a) Each well requires 25  $\mu\text{l}$  of detection antibodies (1x): \_\_\_\_\_ (1)  $\times$  25  $\mu\text{l}$  = \_\_\_\_\_  $\mu\text{l}$  (9)

b) Include a 25% excess to ensure enough volume: \_\_\_\_\_  $\mu\text{l}$  (9)  $\times$  0.25 = \_\_\_\_\_  $\mu\text{l}$  (10)

c) Total volume of 1x detection antibodies: \_\_\_\_\_  $\mu\text{l}$  (9) + \_\_\_\_\_  $\mu\text{l}$  (10) = \_\_\_\_\_  $\mu\text{l}$  (11)

d) Enter the number of diabetes single set (or analytes) tubes that will be multiplexed = \_\_\_\_\_(5)

e) Volume of **20x Detection Antibodies** required from **each diabetes detection antibody tube**:

\_\_\_\_\_  $\mu\text{l}$  (11) /20 = \_\_\_\_\_  $\mu\text{l}$  (12)

f) Total volume of diabetes detection antibody stock: \_\_\_\_\_  $\mu\text{l}$  (12)  $\times$  \_\_\_\_\_(5) = \_\_\_\_\_  $\mu\text{l}$  (13)

g) Volume of **Detection Antibody Diluent** required: \_\_\_\_\_  $\mu\text{l}$  (11) – \_\_\_\_\_  $\mu\text{l}$  (13) = \_\_\_\_\_ $\mu\text{l}$  (14)

## Calculations for Streptavidin-PE

3. Determine the volume of 1x streptavidin PE needed.

a) Each well requires 50  $\mu\text{l}$  of streptavidin PE (1x): \_\_\_\_\_ (1)  $\times$  50  $\mu\text{l}$  = \_\_\_\_\_  $\mu\text{l}$  (15)

b) Include a 25% excess to ensure enough volume: \_\_\_\_\_  $\mu\text{l}$  (15)  $\times$  0.25 = \_\_\_\_\_  $\mu\text{l}$  (16)

c) Total volume of 1x detection antibodies: \_\_\_\_\_  $\mu\text{l}$  (15) + \_\_\_\_\_  $\mu\text{l}$  (16) = \_\_\_\_\_  $\mu\text{l}$  (17)

d) Volume of **100x Streptavidin PE** required: \_\_\_\_\_  $\mu\text{l}$  (17) / 100 = \_\_\_\_\_  $\mu\text{l}$  (18)

e) Volume of **Assay Buffer** required: \_\_\_\_\_  $\mu\text{l}$  (17) \_\_\_\_\_  $\mu\text{l}$  (18) = \_\_\_\_\_  $\mu\text{l}$  (19)

If **multiplexing diabetes (20x) and cytokine (10x) assays**, follow these directions:

Enter the number of wells that will be used in the assay: \_\_\_\_\_ (1)

Enter the number of diabetes tubes (either single set or multiplex) that will be multiplexed: \_\_\_\_\_ (2)

Enter the number of cytokine tubes (either single set or multiplex) that will be multiplexed: \_\_\_\_\_ (3)

## Calculations for Coupled Beads

1. Determine the volume of 1x diabetes and cytokines coupled beads needed.

a) Each well requires 50  $\mu\text{l}$  of coupled beads (1x): \_\_\_\_\_ (1)  $\times$  50  $\mu\text{l}$  = \_\_\_\_\_  $\mu\text{l}$  (4)

b) Include a 20% excess to ensure enough volume: \_\_\_\_\_  $\mu\text{l}$  (4)  $\times$  0.20 = \_\_\_\_\_  $\mu\text{l}$  (5)

c) Total volume of 1x coupled beads: \_\_\_\_\_  $\mu\text{l}$  (4) + \_\_\_\_\_  $\mu\text{l}$  (5) = \_\_\_\_\_  $\mu\text{l}$  (6)

d) **Volume of 20x diabetes coupled beads stock** required from each diabetes tube(s):

\_\_\_\_\_  $\mu\text{l}$  (6) / 20 = \_\_\_\_\_  $\mu\text{l}$  (7)

e) **Volume of 10x cytokines coupled beads stock** required from each cytokines tube(s):

\_\_\_\_\_  $\mu\text{l}$  (6) / 10 = \_\_\_\_\_  $\mu\text{l}$  (8)

f) Total volume of diabetes bead stock required: \_\_\_\_\_  $\mu\text{l}$  (7)  $\times$  \_\_\_\_\_ (2) = \_\_\_\_\_  $\mu\text{l}$  (9)

g) Total volume of cytokine bead stock required: \_\_\_\_\_  $\mu\text{l}$  (8)  $\times$  \_\_\_\_\_ (3) = \_\_\_\_\_  $\mu\text{l}$  (10)

h) Total volume of diabetes and cytokine bead stock required: \_\_\_\_\_  $\mu\text{l}$  (9) + \_\_\_\_\_ (10) = \_\_\_\_\_  $\mu\text{l}$  (11)

i) Volume of **Assay Buffer** required: \_\_\_\_\_  $\mu\text{l}$  (6) \_\_\_\_\_  $\mu\text{l}$  (11) = \_\_\_\_\_  $\mu\text{l}$  (12)

## Calculations for Detection Antibodies

2. Determine the volume of 1x diabetes and cytokines detection antibodies needed.

a) Each well requires 25  $\mu\text{l}$  of detection antibodies (1x): \_\_\_\_\_ (1)  $\times$  25  $\mu\text{l}$  = \_\_\_\_\_  $\mu\text{l}$  (13)

b) Include a 25% excess to ensure enough volume: \_\_\_\_\_  $\mu\text{l}$  (13)  $\times$  0.25 = \_\_\_\_\_  $\mu\text{l}$  (14)

c) Total volume of 1x detection antibodies: \_\_\_\_\_  $\mu\text{l}$  (13) + \_\_\_\_\_  $\mu\text{l}$  (14) = \_\_\_\_\_  $\mu\text{l}$  (15)

d) **Volume of 20x Detection Antibodies** required from each diabetes tube(s):

\_\_\_\_\_  $\mu\text{l}$  (15) / 20 = \_\_\_\_\_  $\mu\text{l}$  (16)

e) **Volume of 10x Detection Antibodies** required from each cytokines tube(s):

$$\text{_____ } \mu\text{l (15)} / 10 = \text{_____ } \mu\text{l (17)}$$

f) Total volume of diabetes detection antibodies stock required: \_\_\_\_\_  $\mu\text{l}$  (16) x \_\_\_\_\_ (2) = \_\_\_\_\_  $\mu\text{l}$  (18)

g) Total volume of cytokine detection antibodies stock required: \_\_\_\_\_  $\mu\text{l}$  (17) x \_\_\_\_\_ (3) = \_\_\_\_\_  $\mu\text{l}$  (19)

h) Total volume of diabetes and cytokine detection antibodies required:

$$\text{_____ } \mu\text{l (18)} + \text{_____ (19)} = \text{_____ } \mu\text{l (20)}$$

i) Volume of **Detection Antibody Diluent** required: \_\_\_\_\_  $\mu\text{l}$  (15) – \_\_\_\_\_  $\mu\text{l}$  (20) = \_\_\_\_\_  $\mu\text{l}$  (21)

## Calculations for Streptavidin-PE

3. Determine the volume of 1x streptavidin PE needed.

d) Each well requires 50  $\mu\text{l}$  of streptavidin PE (1x): \_\_\_\_\_ (1) x 50  $\mu\text{l}$  = \_\_\_\_\_  $\mu\text{l}$  (15)

e) Include a 25% excess to ensure enough volume: \_\_\_\_\_  $\mu\text{l}$  (15) x 0.25 = \_\_\_\_\_  $\mu\text{l}$  (16)

f) Total volume of 1x detection antibodies: \_\_\_\_\_  $\mu\text{l}$  (15) + \_\_\_\_\_  $\mu\text{l}$  (16) = \_\_\_\_\_  $\mu\text{l}$  (17)

d) Volume of **100x Streptavidin PE** required: \_\_\_\_\_  $\mu\text{l}$  (17) / 100 = \_\_\_\_\_  $\mu\text{l}$  (18)

e) Volume of **Assay Buffer** required: \_\_\_\_\_  $\mu\text{l}$  (17) \_\_\_\_\_  $\mu\text{l}$  (18) = \_\_\_\_\_  $\mu\text{l}$  (19)