

MULTISCREEN™ Calcium 1.0 No-Wash Assay Kit Protocol

INTRODUCTION

Multiscreen™ Calcium 1.0 No Wash Assay Kits provide homogeneous fluorescence-based assays for intracellular calcium mobilization detection in 96-well, 384-well and higher multiplexity. Pre-loaded with Calcium 1.0 is taken up by the cells where the lipophilic blocking groups of Calcium 1.0 are cleaved by esterases, resulting in a negatively charged fluorescent dye that stays inside cells. The fluorescence signal is greatly enhanced after binding to intracellular calcium released by cell-stimulation. Long wavelength, high sensitivity, and >100 times fluorescence enhancement make Calcium 1.0 an ideal indicator for intracellular calcium release assays measuring GPCRs, calcium channels and receptor tyrosine kinase signaling. Multiscreen™ Calcium 1.0 No Wash Assay Kits are validated by Multiscreen™ stable cell lines.

Kit Components <i>(Warning: Do not add additional probenecid)</i>	Catalog Numbers			Storage	Instruments
	MSCA01-1 (1K tests, 384-well)	MSCA01-10 (10K tests, 384-well)	MSCA01-100 (100K tests, 384-well)		
Component A: Calcium 1.0 + probenecid	1 vial, lyophilized	1 vial, lyophilized	10 vials, lyophilized	-20°C and avoid light	FLIPR, FDSS NOVOStar FlexStation ViewLux ArrayScan
Component B: 10X Pluronic® F127 Plus	1 bottle (1 mL/bottle)	10 bottles (1 mL/bottle)	10 bottles (10 mL/bottle)		
Component C: HHBS (HANKS + 20 mM HEPES)	1 bottle (9 mL)	1 bottle (100 mL)	Not included		

CALCIUM ASSAY PROTOCOL

1. Prepare cells *(Evaluate each cell line to determine optimal cell density and other conditions.)*

- 1.1 **Adherent:** Plate cells overnight in growth media with 10% FBS at 3,000 to 9,000 cells/well/40 µL for a 384-well black poly-D lysine coated clear bottom plate.
- 1.2 **Suspension:** Centrifuge the cells from the culture medium and then suspend the cell pellet in the appropriate amount of Component C 10,000 to 20,000 cells/well/30 µL for a 384-well black poly-D lysine coated clear bottom plate. Allow cells to settle for 10 minutes at RT. Centrifuge the plate at 200x g for 5 minutes with brake off prior to the experiments.

2. Prepare Calcium 1.0 dye-loading solution *(Mix well by gentle-vortexing after each step 2.2 - 2.4.)*

- 2.1 Thaw Component A, Component B and Component C at room temperature.
- 2.2 **Make Calcium 1.0 Stock Solution:** Add 20 µL of DMSO into the vial of Component A in the MSCA01-1 kit or 200 µL to each vial of Component A in the MSCA01-10 and MSCA01-100 kits. *Mix well.*
- 2.3 **Make 1X Assay Buffer:**
 - For Catalog MSCA01-1 and MSCA01-10, add 9 mL of Component C into 1 mL Component B. *Mix well*
 - For Catalog MSCA01-100, add entire bottle 10mL of Component B into 90 mL of Component C. Prepare in a separate polypropylene bottle (not included in the kit). *Mix well.*
- 2.4 **Make Calcium 1.0 dye-loading solution:**
 - For Catalog MSCA01-1 and MSCA01-10, add 20 µL of Calcium 1.0 Stock Solution (2.2) to 10 mL of 1X Assay Buffer (2.3). *Mix well.*
 - For Catalog MSCA01-100, add 200 µL of Calcium 1.0 Stock Solution (2.2) to 100 mL of 1X Assay Buffer (2.3). *Mix well.*
 - Aliquot into 10 mL light blocking opaque bottles. Store at -20°C, protect from light, and avoid repeated freeze thaw cycles.
 - The final solution is stable for 2 hours at room temperature.

3. Run Calcium assay:

- 3.1 For adherent cell assay, replace cell growth media with 30 µL/384-well HHBS buffer to minimize background fluorescence and compound interference by serum.
- 3.2 Add Calcium 1.0 dye-loading solution at 10 µL/384-well.
- 3.3 Incubate the dye-loading plate in a cell incubator for 1 hour at 37°C.
- 3.4 Prepare compound plate with Component C
- 3.5 Run Calcium flux assay by monitoring the fluorescence intensity at Ex/Em = 490/ 525 nm.

Note: Prepare instrument by adjusting the signal intensity to 10%-15% of the maximum instrument counts. For example, the maximum for FLIPR-384 is 65,000, so the settings should be adjusted to 7,000 to 10,000.

