

MULTISCREEN™ TR-FRET cAMP 1.0 No-Wash Assay Kit

INTRODUCTION

Multiscreen™ TR-FRET cAMP 1.0 No Wash Assay Kits provide a homogenous TR-FRET assay method for adenylyl cyclase activity detection. α -cAMP antibody (Ab) is labeled with MultiScreen™Eu while cAMP is labeled with MultiScreen™650. In the absence of cAMP, MultiScreen™650-cAMP is bound to MultiScreen™Eu- α -cAMP-Ab to give a strong TR-FRET Emission at 655 nm. Free cAMP in the test sample competes for binding to the MultiScreen™ Eu- α -cAMP-Ab, reducing TR-FRET signal from MultiScreen™650-cAMP binding. The MultiScreen™650 labeled cAMP only has fluorescence lifetime of nanosecond while MultiScreen™Eu-labeled α -cAMP antibody has much longer fluorescence lifetime value due to the TR-FRET. The magnitude of TR-FRET is proportional to the concentration of cAMP in a sample. The assay can be performed in 96- or 384-well microtiter-plate format and adapted to automation.

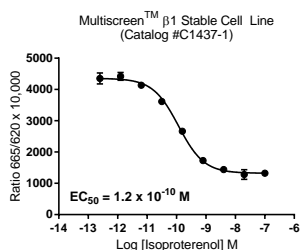
Kit Components	Catalog Numbers			Storage	Instrument Platform
	MSCM01-1 (0.5K tests, 384-well)	MSCM01-10 (5K tests, 384-well)	MSCM01-100 (25K tests, 384-well)		
Component A: MultiScreen™Eu- α -cAMP-Ab	1vial	1 vial 50x	5 vials 50x	-4 °C and avoid light	TR-FRET microplate readers
Component B: MultiScreen™650-cAMP	1vial	1 vial 50x	5 vials 50x		
Component C: cAMP Standard	1 vial (20 μ L 1mM)	1 vial (33 μ g)	1 vial (33 μ g)		
Component D: Cell Lysis Buffer	10 mL	100 mL	100 mL 5 bottles		
Component E: Diluent	10 mL	100 mL	100 5 bottles		
Reagents NOT included in the kit	Assay Buffer: Hank's Balanced Salt Solution with 20mM HEPES, pH 7.4; When necessary: IBMX (3-Isobutyl-1-methylxanthine); Forskolin (Adenylyl cyclase activator)				

PREPARE WORKING SOLUTIONS (Mix well by gentle-vortexing or pipette mixing after each step):

- cAMP standard:** Add 100 μ L Component E to Component C to make 1mM stock solution for MSCM01-10 and MSCM01-100 kits only and gently vortex to mix: Add 1 μ L 1mM stock solution into 99 μ L Component E or cell culture media to make 10 μ M standard followed 4-fold serial dilution to make 10000, 2500, 625, 156.3, 39.1, 9.8, 2.4, 0.61, 0.15, 0.038 nM final concentrations. Mix gently with a pipette after each dilution. Add 10 μ L or 20 μ L of serial diluted cAMP standard per 384-well in microtiter assay plate after last the incubation period from step 5.
- MultiScreen™Eu- α -cAMP-Ab working solution:** Add 50 μ L Component A to 2.5mL Component D (scale down based on need). Prepare right before use. Store at 4°C.
- MultiScreen™650-cAMP working solution:** Add 50 μ L Component B to 2.5mL Component D (scale down based on need). Prepare right before use. Store at 4°C.

cAMP ASSAY PROTOCOL (384-well format)

- Prepare cells** (Evaluate each cell line to determine optimal cell density and other conditions.)
Adherent: Plate cells overnight in growth media with 10% FBS at 3k-9k cells/40 μ L/384-well in Poly-D-Lysine coated white opaque bottom plate. Remove growth media carefully before compound treatment.
Suspension: Centrifuge the cells from the culture media and then suspend the cell pellet in the appropriate amount of Assay Buffer at 3k-12k cells/5 μ L/ 384-well in small volume, white, opaque bottom plate.
- Compound Treatment** (The incubation time and temperature can be optimized for each receptor): Add 15 μ L of test compounds to adherent cells or 5 μ L to suspension cells per well and incubate for 20 minutes at 37°C.
- Termination (30 μ L or 20 μ L final volume):** Add 7.5 μ L MultiScreen™650-cAMP working solution and 7.5 μ L MultiScreen™Eu- α -cAMP-Ab working solution sequentially to adherent cells per well or 5 μ L of each on sequentially to suspension cells per well. Incubate 30 minutes at room temperature in the dark. Read fluorescence emission on a TR-FRET compatible reader at 665 nm and 620 nm.



Sample Data: cAMP Dose Response Curve in Multiscreen™ β 1-HEK293T stable cell line (Multispan C1437-1). Measured with MultiScreen™ TR-FRET cAMP 1.0 No Wash Assay Kit. β 1-HEK293T cells were suspended in 5 μ L HBSS assay buffer per well in a 384-well solid white plate. The cells were treated with Isoproterenol and 1mM IBMX for 20 minutes at 37°C. The reaction was terminated with 5 μ L of MultiScreen™Eu- α -cAMP-Ab working solution and 5 μ L MultiScreen™650-cAMP working solution followed by 30-minute incubation at RT. Relative Fluorescence Units were measured at 665nm and 620nm. Ratio is calculated as the F_{665nm} / F_{620nm} ratio and expressed in Delta F% $R = F_{665nm} / F_{620nm}$; Delta F% = 100% x $(R_{sample} - R_{neg}) / R_{neg}$. Standard curve was drawn by plotting Delta F% versus cAMP concentration.