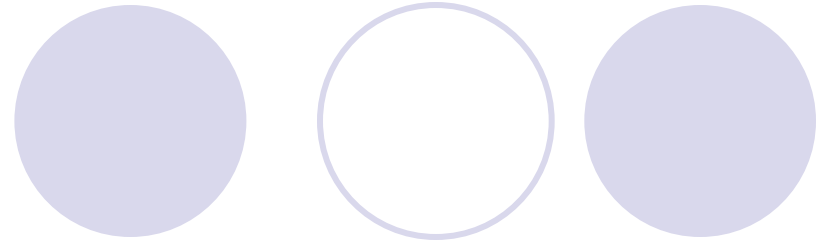
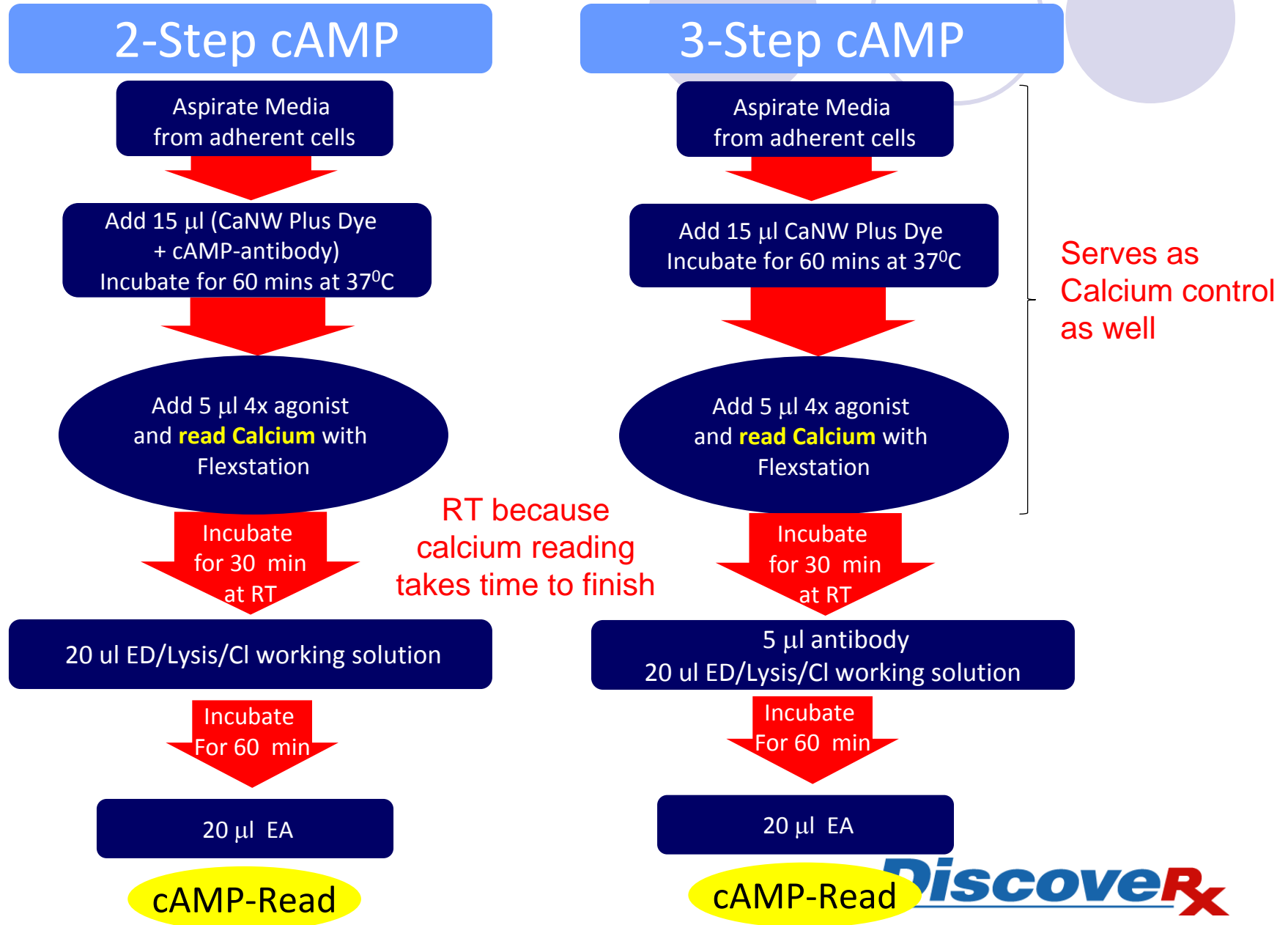


## Purpose of experiment

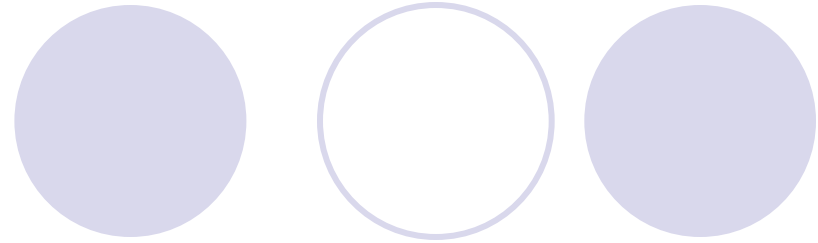


- Feasibility of multiplexing calcium and cAMP read out. To do this we used:
  - Human Prokineticin receptor 1 CHO (hPROKR1) cAMP Hunter cell line (95-0090C2) (Gs coupled)
  - Calcium NO WASH PLUS assay kit (90-0091)
  - cAMP XS+ kit (90-0075)
- Concentration response curve of reference agonist hu EG-VEGF (Peprotech cat# 100-44, lot #0403244)
- Both calcium and cAMP assays were run as a control in parallel to compare results with multiplexing data

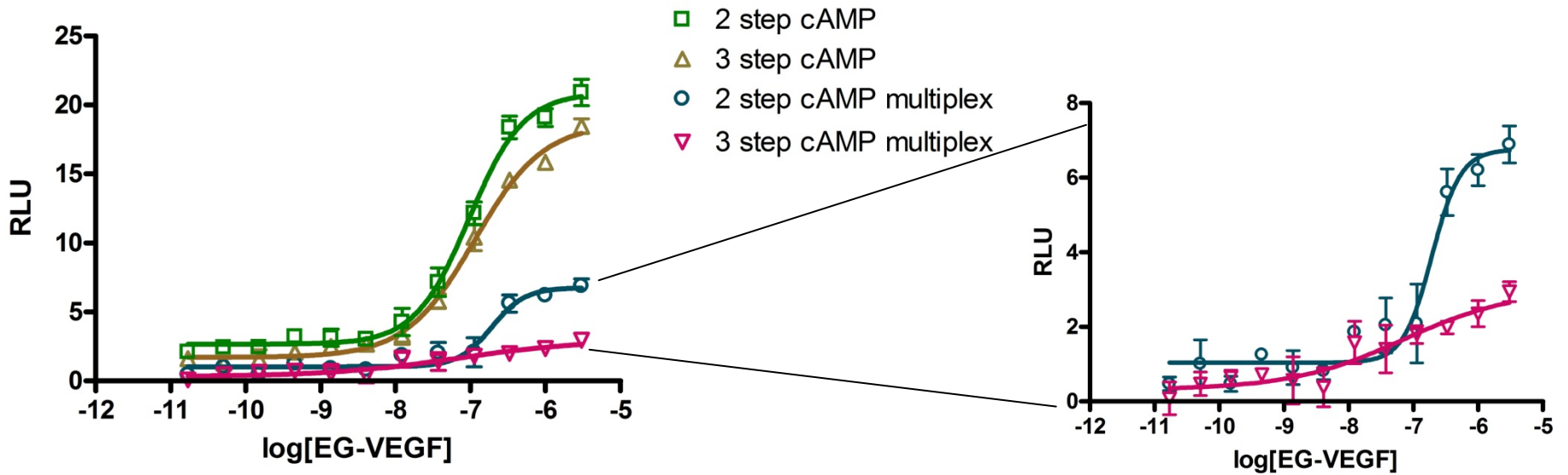
# Experimental set up for MULTIPLEXING



# cAMP



cAMP data on PROKR1



	2 step cAMP	3 step cAMP	2 step cAMP multiplex	3 step cAMP multiplex
BOTTOM	2.655	1.715	1.041	0.3170
TOP	20.89	18.67	6.752	3.000
LOGEC50	-7.018	-6.946	-6.717	-7.159
HILLSLOPE	1.204	0.9479	1.964	0.5011
EC50	9.5981e-008	1.1326e-007	1.9200e-007	6.9308e-008
<b>S:B</b>	<b>8</b>	<b>11</b>	<b>6</b>	<b>~ 3</b>