

Comparison of cAMP Assay Technologies for High Throughput Screening

Patricia Kasila and Harry Harney

PerkinElmer Life Sciences, 549 Albany Street, Boston, MA 02118

1

Abstract

Homogeneous cAMP assays have been developed to allow for the direct measurement of receptor mediated adenylyl cyclase activation/inhibition in G-protein coupled receptor systems. PerkinElmer has developed three homogenous cAMP assays using different technologies. These assays can all be easily modified to accommodate many robotic systems and situations in fully homogeneous assays. The platforms described here include scintillation proximity technology using the Adenylyl Cyclase Activation FlashPlate kit (Cat # SMP701), fluorescence polarization using [FP]² cAMP kit (Cat # FPA202), and luminescent proximity technology using the AlphaScreen cAMP assay (Cat # 6760600C). Studies were performed to compare sensitivity, precision, and ease of use of each assay. Requirements for technology attributes vary greatly depending on the individual user. The comparative data shown here will highlight advantages of each technology.

2

Introduction

cAMP Technology Comparison

- ▶ Adenylyl Cyclase Activation FlashPlate Assay
- ▶ [FP]² cAMP Assay
- ▶ AlphaScreen cAMP Assay

Comparisons

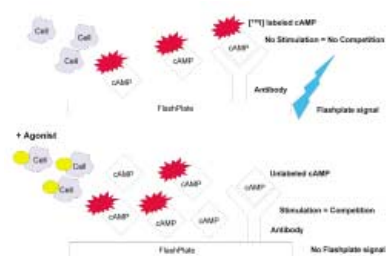
Using whole cells with overexpressed $\beta 2$ Adrenergic receptors

- ▶ Standard Curve Dynamic Range
- ▶ Agonist Dose Response
- ▶ Forskolin Dose Response
- ▶ Antagonist Dose Response

3

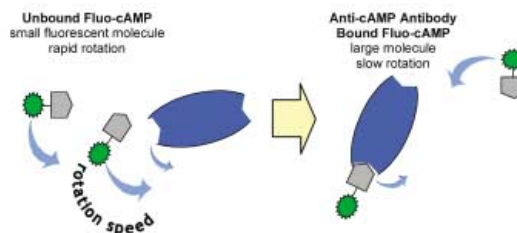
Basic Principle

Adenylyl Cyclase Activation FlashPlate Assay



- ▶ agonist-stimulated cAMP displaces [¹²⁵I] labeled cAMP from Antibody

[FP]² Fluorescence Polarization cAMP Assay

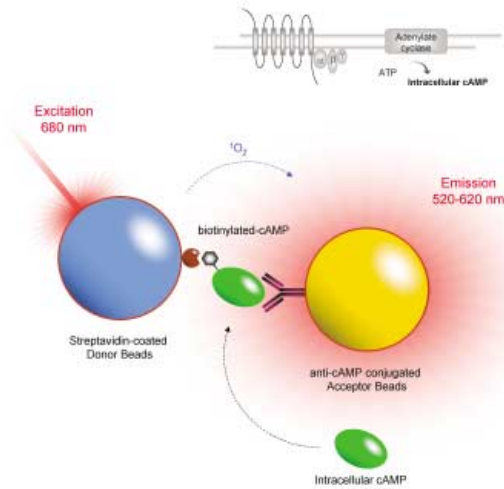


- ▶ agonist-stimulated cAMP displaces Fluo-cAMP from Antibody produces lower polarization signal

4

Basic Principle (cont'd)

AlphaScreen cAMP



- › agonist-stimulated cAMP displaces biotinylated cAMP from anti-cAMP immobilized on acceptor beads.

5

Kit Components

Adenylyl Cyclase Activation FlashPlate Assay (PerkinElmer Catalog # SMP701)

- › cAMP Standard
- › Stimulation Buffer
- › [¹²⁵I]cAMP Tracer
- › Detection Buffer
- › Adenylyl Cyclase FlashPlates

[FP]² Fluorescence Polarization cAMP Assay (PerkinElmer Catalog # FPA202)

- › cAMP Standard
- › Stimulation Buffer
- › cAMP Antibody
- › Fluo-cAMP Tracer
- › Detection Buffer
- › Microtiter plates

AlphaScreen cAMP Assay (PerkinElmer Catalog # 6760600C)

- › Biotinylated cAMP
- › Streptavidin Donor Beads
- › Anti-cAMP Acceptor Beads

6

Assay Protocols

Adenylyl Cyclase Activation FlashPlate Assay

- 1- 10 μ L of cAMP standards or cells diluted in Stimulation Buffer
- 2- 15 μ L PBS (standards) or forskolin or agonist/antagonist (cells)
- 3- Incubate for 30 – 60 min. at room temperature
- 4- 25 μ L of Detection Mix containing Detection Buffer and [¹²⁵I]cAMP
- 5- Incubate for 2 hours at room temperature
- 6- Read on a MicroBeta[®] or TopCount NXT[®] Microplate Scintillation Counter

[FP]² cAMP

- 1- 10 μ L of cAMP standards or cells diluted in Stimulation Buffer
- 2- 10 μ L PBS (standards) or forskolin or agonist/antagonist (cells)
- 3- Incubate for 30 – 60 min. at room temperature
- 4- 10 μ L of Detection Mix containing Detection Buffer and Fluo-cAMP (5 nM final)
- 5- 10 μ L of Anti-cAMP diluted in Detection Buffer
- 6- Incubate for 60 minutes at room temperature
- 7- Read on a Fluorescence Polarization Reader (EnVision[™], ViewLux[™], Victor²V[™] or Fusion[™])

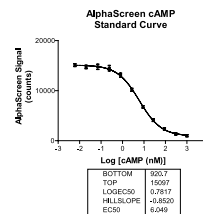
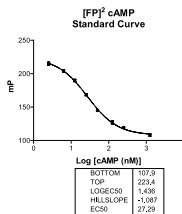
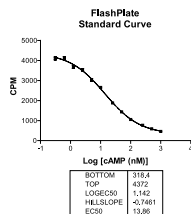
AlphaScreen cAMP

- 1- 5 μ L cells/anti-cAMP Acceptor Beads or anti-cAMP Acceptor Beads only (standards)
- 2- 5 μ L forskolin or agonist/antagonist or cAMP standards
- 3- Incubate for 30 – 60 min. at room temperature
- 4- 15 μ L of biotinylated-cAMP/Streptavidin Donor Beads Detection Mix
- 5- Incubate for at least 60 min. at room temperature
- 6- Read on a Fusion[™]- α or AlphaQuest[™]-HTS microplate analyzer

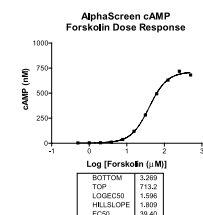
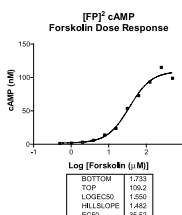
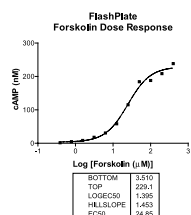
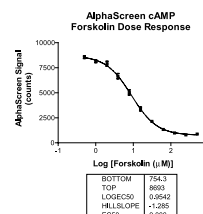
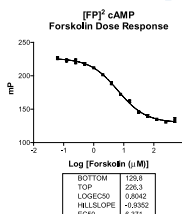
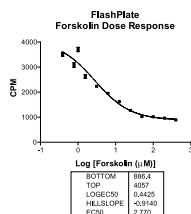
7

Results

cAMP Standard Curves

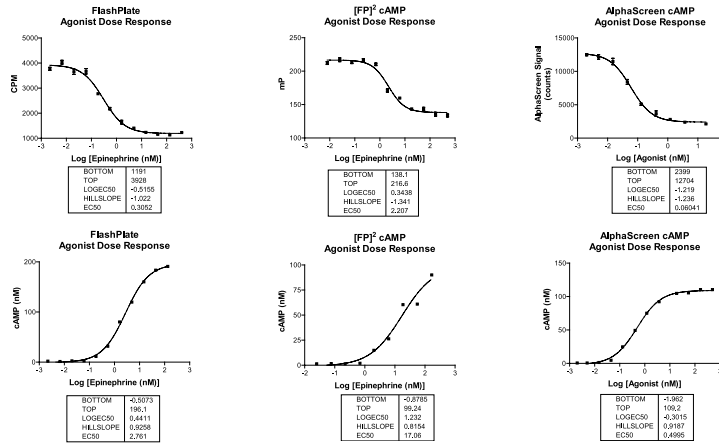


Forskolin Dose Response

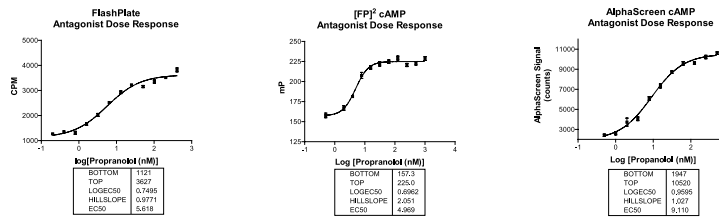


Results (cont'd)

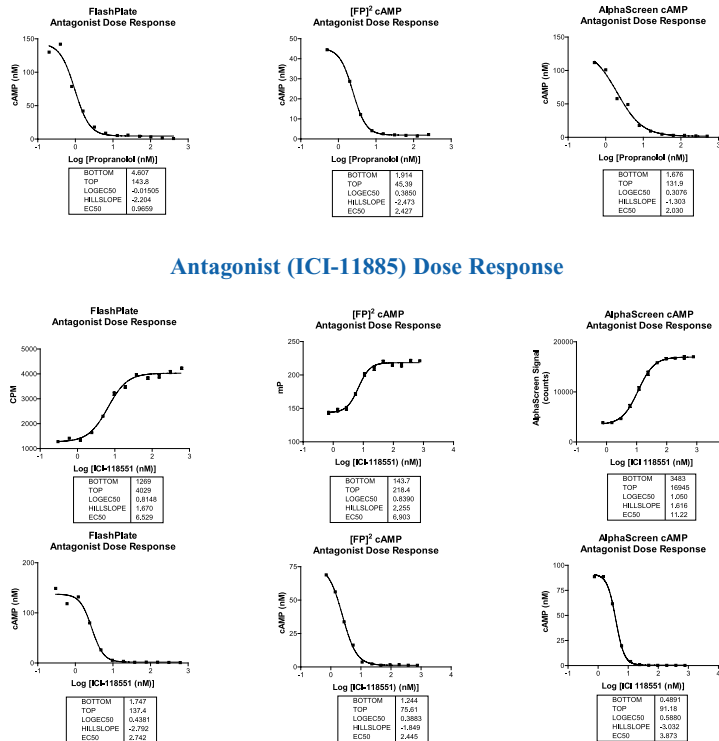
Agonist Dose Response



Antagonist (Propranolol) Dose Response



Antagonist (ICI-11885) Dose Response



8

Results (cont'd)

EC₅₀ and Z' Comparison (Raw Data)

	Standard Curve		Forskolin Dose Response		Agonist Dose Response		Antagonist Dose Response (Propranolol)		Antagonist Dose Response (ICI-11885)	
	EC ₅₀	Z'	EC ₅₀	Z'	EC ₅₀	Z'	EC ₅₀	Z'	EC ₅₀	Z'
FlashPlate	13.9	0.98	2.8	0.97	0.31	0.94	5.6	0.84	6.5	0.92
[FP] ²	27.3	0.87	6.4	0.74	2.2	0.80	5.0	0.76	6.9	0.87
AlphaScreen	6.0	0.77	9.0	0.70	0.06	0.89	9.1	0.86	11.2	0.84

EC₅₀ of Interpolated Results

	Forskolin Dose Response	Agonist Dose Response	Antagonist Dose Response (Propranolol)	Antagonist Dose Response (ICI-11885)
FlashPlate [cAMP(nM)]	24.8	2.8	0.97	2.7
[FP] ² [cAMP(nM)]	35.5	17.1	2.4	2.4
AlphaScreen [cAMP(nM)]	39.4	0.50	2.0	3.9

9

Conclusions

The cAMP technologies presented here are homogeneous assays that are highly suited for High Throughput Screening. The Adenylyl Cyclase FlashPlate Assay has been used in many HTS screening labs and has been found to be the “gold standard” for cAMP detection. [FP]² cAMP fluorescence polarization assay is a non-rad homogenous platform that is easy to use, and gives the flexibility to format the protocol depending on the needs of the user. This assay does not have the dynamic range or sensitivity of the other technologies but can be used successfully for many receptors. AlphaScreen cAMP has the widest dynamic range, excellent precision characteristics and is a valid non-rad alternative to the “gold standard” Adenylyl Cyclase FlashPlate assay.



Worldwide Headquarters: PerkinElmer Life Sciences, Inc., 549 Albany Street, Boston, MA 02118-2512 USA (800) 551-2121

European Headquarters: PerkinElmer Life Sciences, Inc., Imperiastraat 8, BE-1930 Zaventem Belgium

Technical Support in Europe: techsupport.europe@perkinelmer.com in US and Rest of World: techsupport@perkinelmer.com

Belgium: Tel: 0800 94 540 • **France:** Tel: 0800 90 77 62 • **Netherlands:** Tel: 0800 02 23 042 • **Germany:** Tel: 0800 1 81 00 32 • **United Kingdom:** Tel: 0800 89 60 46
Switzerland: Tel: 0800 55 50 27 • **Italy:** Tel: 800 79 03 10 • **Sweden:** Tel: 020 79 07 35 • **Norway:** Tel: 800 11 947 • **Denmark:** Tel: 80 88 3477 • **Spain:** Tel: 900 973 255