

High Throughput Tag-Lite® Cell-based Functional and Surface Binding Assays on the SpectraMax® Paradigm Plate Reader Platform

Oksana Sirenko¹, Marie-Laure Lebreton², Delphine Jaga², Michael Katzlinger¹, Laurence Monnet¹, Caroline Cardonnel¹, H. Roger Tang¹, Cathy Olson¹, Francois Degorce², Evan F. Cromwell¹

¹Molecular Devices, Inc. Sunnyvale, CA

²Cisbio Bioassays, Codolet, France

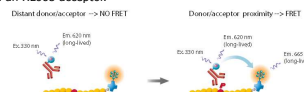
Introduction

G-protein coupled receptors (GPCR) are the largest class of cell-surface receptors and are targets for almost 40% of existing drugs. Lead discovery, testing the efficacy of prospective drugs (in the area of cardiovascular diseases and other fields), and understanding of mechanism of action of drug candidates requires assays that can measure the binding of ligands to the receptors, receptor oligomerization, and/or internalization. Accordingly, there is a real need for robust and sensitive assays of this type that are suitable for high throughput screening. The Tag-Lite® cellular screening platform was designed to increase the flexibility of cell-surface receptor research. This platform, which combines HTRF® and SNAP-tag® technologies, is ideal for primary and secondary screening and can be applied to a variety of assay formats for pharmacological characterization and development of therapeutic antibodies. Here we show results from use of this assay platform on the only user upgradable plate reader – the SpectraMax® Paradigm system, and the widely used M5e plate reader.

Method

Principles of TR-FRET measurements

HTRF® technology (Homogeneous Time-Resolved Fluorescence) is a TR-FRET based read out that uses the principles of both TRF and FRET. The HTRF donor fluorophore is either Europium cryptate (Eu3+ cryptate) or Lumi4®-Tb (Tb2+ cryptate), fruit of a recent collaboration with Lumipore Inc. Various acceptor molecules can be used which are either Red or Green emitters. When the two fluorophores are brought together by a biomolecular interaction, a portion of the energy captured by the Cryptate during excitation is released through the acceptor. The cartoon below shows the basic principles of TR-FRET using an Eu3+ cryptate donor and an XL665 acceptor.



SpectraMax® Paradigm and M5e Plate Reader Platforms



- SpectraMax® Paradigm System**
- User upgradeable high throughput multi-mode reader w/dual PMTs
 - High sensitivity for all applications
 - Accepts all standard microplates up to 1536 wells
 - The Paradigm HTRF cartridge was used for Validation, pAKT, pERK, and Tag-Lite binding assays
 - Ex 330nm, Em 616nm & 665nm
 - The Paradigm Tb-Green cartridge was used for the cAMP assay
 - Ex 330nm, Em 490nm & 520nm
 - Read time for 384 well plate: 2.3 min

- SpectraMax® M5e System**
- Five modes of detection for wide range of applications
 - The standard for UV/Vis absorbance
 - SoftMax® Pro industry leading, all-in-one plate reader software
 - Setup for Validation, pAKT, pERK, and Tag-Lite binding assays:
 - Ex 330nm, Em1 616nm, Em2 665nm
 - Setup for cAMP assay:
 - Ex 330nm, Em1 490nm, Em2 520nm

Results and Discussion

cAMP Detection Assay

cAMP assay kits allow direct quantitative determination of cyclic AMP with either suspended or adherent cells using HTRF reagents. The method employs a competitive immunoassay between native cAMP produced by cells and the cAMP labeled with the dye d2. The tracer binding is visualized by a Mab anti-cAMP labeled with the dye d2. The capability of the assay on the SpectraMax Paradigm and M5e readers was evaluated by titration curves shown below.

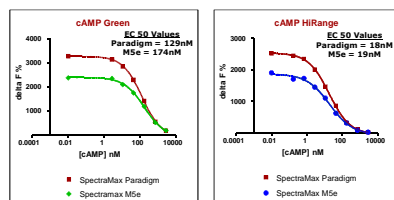
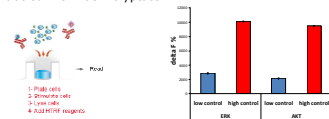


Figure 1. Cyclic AMP titration curves were evaluated using SpectraMax Paradigm and SpectraMax M5e instruments. Both instruments demonstrated good performance for the assay. Paradigm: W = 21.4, Z' = 0.91. M5e: W = 15.5, Z' = 0.97

HTRF Cellular Kinase Assays for phospho-AKT and phospho-ERK

Cellular erk and HTRF phospho-Akt (Ser473) assays allow detection of activated Erk1/2 and Akt directly in whole cells. Upon receptor activation, the kinases are activated, and upon cell lysis phosphorylated kinases can be detected using the kit reagents. The assays are based on a sandwich immunoassay involving anti-kinase antibody labeled with d2, and an anti-phospho-kinase antibody labeled with Eu3+ cryptate.



SpectraMax Paradigm		SpectraMax M5e	
ERK	AKT	ERK	AKT
low control	high control	low control	high control
3872	10462	2104	8626
276	1%	25	4%
3.5	4.4	3.3	3.6

Figure 2. Stimulated and un-stimulated cell lysates provided as assay internal controls to check the quality of the results obtained. The window between high and low controls, shown in the bottom line of the table, should be greater than 2.

Validation of Plate Readers for Cisbio Assays:

The HTRF ratio is calculated by $\text{Ratio} = \frac{A_{620nm}}{B_{620nm}} \times 10^4$. ΔF is used to improve assay reproducibility $\Delta F = \frac{\text{Ratio}_{low} - \text{Ratio}_{high}}{\text{Ratio}_{high}} \%$. W is used to define the Assay Window $W = \frac{\text{Ratio}_{high}}{\text{Ratio}_{low}}$.

Standard d	Parameter	Requirement	Paradigm	M5e
Low Calibrator	CV Ratio	<10%	5%	7%
	ΔF	>15%	28%	21%
High Calibrator	ΔF	>600%	910%	664%
S/B		>40	96	50

Table 1: Results for SpectraMax Paradigm and M5e plate reader certification tests in 384 well white plates. Paradigm system has one of the highest S/B ratios for HTRF among multimode readers.

Tag-Lite® Live Cell GPCR Binding Assays

The Tag-Lite platform allows one to efficiently label a protein of interest on a targeted site with HTRF dyes. Cisbio Bioassays offers plasmids encoding TAGS and protein of interest, or frozen cells already transfected with the constructs. The constructs lead to the expression of the tagged protein that can be labeled with Terbium Cryptate, while the receptor ligand (agonist or antagonist) is conjugated with acceptor. The Tag-Lite platform is ideal for a wide range of applications, such as mechanistics and receptor dimerization, ligand binding assays, and second messenger assessment.

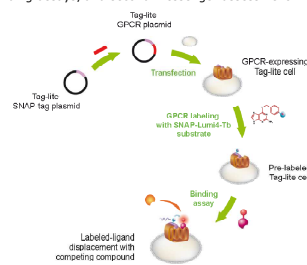


Figure 3. Depiction of the Tag-Lite cell surface binding assay protocol.

Optimization of instrumental settings for SpectraMax Paradigm and SpectraMax M5e

Delay time and Integration times were modified during optimization. Greater assay window was achieved when shorter delay and integration times were used, for both instruments.

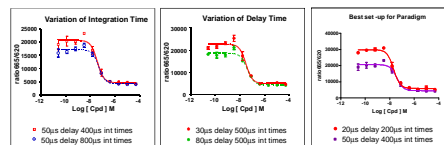


Figure 4. Characteristic titration curves from the assay optimization experiments. Comparison was made of different integration times (Left) and delay times (Middle). The optimized response curve (20µs delay, 200µs int) is shown in the plot on the Right.

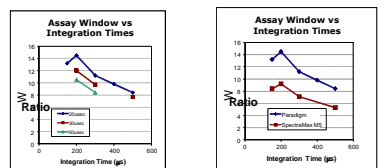


Figure 5. Plots of the assay window value W for various integration and delay times. Left: Assay optimization results for the Paradigm system. The optimum settings were found to be 20µs delay, 200µs int. Right: Comparison between SpectraMax Paradigm and M5e plate readers. The systems showed similar optimization behavior.

Dopamine D3 Binding Assay:

Dopamine receptors are a class of G protein-coupled receptors that are prominent in the vertebrate central nervous system. Dopamine receptors are implicated in many neurological processes, including pleasure, cognition, memory, and fine motor control. Abnormal dopamine receptor signaling is implicated in neuropsychiatric disorders, thus dopamine receptors are common neurologic drug targets. Antipsychotic drugs are often dopamine receptor antagonists, while psychostimulants are typically indirect agonists of dopamine receptors. A D3 cell-based Tag-Lite binding assay allows testing of receptor-selective agonists and antagonists and possibly development of novel antipsychotic drugs.

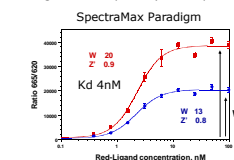


Figure 6. Binding assay: HEK293 cells expressing D3 receptor were labeled with Tb Cryptate, then plated and incubated with donor-conjugated ligand for 90min. Different delay and integration times for both instruments were assessed. Shorter integration time resulted in greater W and Z' values.

We have evaluated performance of the Dopamine D3 Tag-Lite binding assay for the SpectraMax Paradigm and SpectraMax M5e plate readers using optimized settings. Results from a competitive inhibition assay on the two systems are shown below. The Paradigm reader shows superior results and offers unmatched flexibility. Users can purchase assay-optimized cartridges and upgrade the system at any time to enable capabilities for future applications

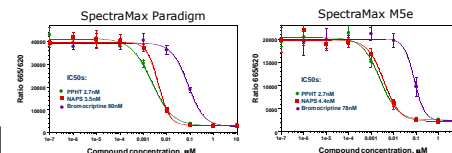


Figure 7. Competitive inhibition assay: several known dopamine receptor agonists and antagonists were tested and IC50s determined. Cells were incubated in the presence of 6nM of donor-conjugated ligand as well as one of the receptor agonists PPH1 and bromocriptine or antagonist NAGS. Similar IC50s were determined in the assay by both plate readers.

Summary

- While both the Paradigm and M5e plate readers demonstrated excellent performance for all tested assays, the Paradigm plate reader demonstrated better sensitivity and superior performance for all of these assays.
- Excellent performance of the SpectraMax Paradigm and M5e plate readers was demonstrated for several Cisbio assays: cAMP detection, pERK and pAKT kinase assays, and Tag-Lite Dopamine D3 receptor ligand binding assay.
- Optimization of instrumental settings was done on both instruments for the Tag-Lite assays in 384 multi-well plate format. Performance was evaluated by both assay window and Z-prime values.
- Shortening both delay and integration times resulted in better instrument performance for the SpectraMax Paradigm plate reader.
- The combination of the SpectraMax Paradigm plate reader with the Tag-Lite functional assay is shown to be a powerful assay platform for cell surface binding and functional assays and well suited for high throughput screening.



FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES. Molecular Devices, the Molecular Devices logo, and all other trademarks, unless otherwise indicated, are the property of Molecular Devices, Inc. ©2011 Molecular Devices, Inc. HTRF and Tag-Lite are registered trademarks of Cisbio Bioassays. SNAP tag is a registered trademark of New England Biolabs.