

# HTS instrument discovers low affinity inhibitors of the Inositol Phosphate (IP) signaling pathway

EJ Dell<sup>1</sup>, JL Tardieu<sup>2</sup>, and F Degorce<sup>2</sup>  
<sup>1</sup>BMG LABTECH GmbH <sup>2</sup>Cisbio Bioassays

- Several low affinity drug 'hits' were only found by the PMT-based reader and not the CCD-based reader
- PHERAstar<sup>®</sup> FS performs more than twice as fast as the CCD-based reader
- PHERAstar<sup>®</sup> FS shows superior assay quality parameters such as Z', DeltaF% and assay window

## Introduction

G-protein coupled receptors (GPCRs) are transmembrane proteins which play a key role in the signal transduction of extracellular stimuli. GPCRs are associated with a complex assembly of intracellular proteins regulating a large variety of downstream effectors. The production of inositol 1,4,5-triphosphate (IP<sub>3</sub>) is one such second messenger, which is produced in response to the activation of Gq-coupled receptors. However, IP<sub>3</sub>'s very short half-life make its assessment too challenging for drug screening assays and the monitoring of calcium release, triggered by IP<sub>3</sub>, has been extensively used as a downstream readout of this signaling pathway. An alternate way to monitor IP<sub>3</sub> is to measure the accumulation of inositol monophosphate (IP<sub>1</sub>), which is a downstream metabolite of IP<sub>3</sub>.

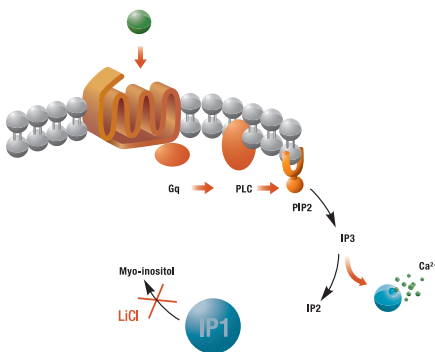


Fig. 1: Biosynthesis of IP<sub>1</sub>.

Monoclonal antibodies raised against IP<sub>1</sub> lead to the optimization of a homogeneous time-resolved fluorescence assay [HTRF<sup>®</sup>], taking advantage on the fact that IP<sub>1</sub> is stable and accumulates in cells. The HTRF<sup>®</sup> IP-One assay has been compared to existing methods and has been shown to lead to similar compound potency data. Moreover, the end point accumulation of IP<sub>1</sub> allows for the discrimination of slow acting compounds that remain unseen by calcium sensing. The HTRF<sup>®</sup> IP-One assay also allows for the characterization of inverse agonists by the quantification of constitutively active GPCRs, which is impossible via measurement of calcium release. Lastly, HTRF<sup>®</sup> IP-One detection confers superior assay robustness and much lower false positive rates compared to calcium detection. The IP-One assay developed by Cisbio Bioassays uses their proprietary HTRF<sup>®</sup> technology (Figure 2).

## Assay Principle

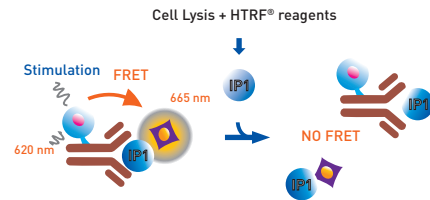


Fig. 2: IP-One HTRF<sup>®</sup> Assay Principle.

The assay uses a monoclonal antibody that specifically recognizes IP<sub>1</sub> and it is based on a competition format in which the intracellular accumulation of IP<sub>1</sub> inhibits the fluorescence resonance energy transfer (FRET) signal between the HTRF<sup>®</sup> donor and acceptor. An IP<sub>1</sub> calibration curve can estimate the IP<sub>1</sub> concentration accumulated in cells as a function of the compound concentration.

## Materials & Methods

- Next generation PMT based HTS microplate reader, PHERAstar<sup>®</sup> FS
- HTS CCD-based microplate imager from a different vendor

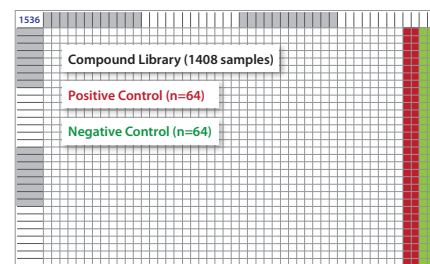


Fig. 3: 1536-well microplate layout.

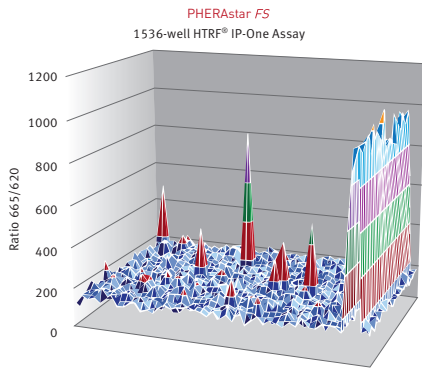
Black Greiner 1536-well microplates were used with an assay volume of 5  $\mu$ L. As shown in Figure 3, 1408 different compounds were pipetted into the first 44 columns of a microplate, with positive (POS, n=64) and negative (NEG, n=64) controls in the last 4 columns. The PHERAstar<sup>®</sup> FS is equipped with a high power pulsed nitrogen laser emitting at 337 nm, as well as a dedicated Simultaneous Dual Emission (SDE) direct photon counting time-resolved fluorescence mode. When exciting the terbium (Tb) donor molecule, the



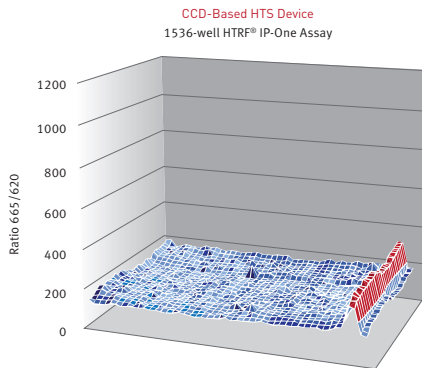
laser is superior to a broadband xenon flash lamp. The laser's energy emission takes advantage of the higher molecular extinction coefficient of the terbium cryptate peaking around 337 nm, compared to europium [Eu].

## Results & Discussion

Results were evaluated by the HTRF<sup>®</sup> ratio of the two emission wavelengths (Em 665 nm / Em 620 nm) and hits are shown as peaks in a surface graph representing the entire 1536 well microplate.



**Fig. 4:** HTRF<sup>®</sup> ratios obtained for the IP-One assay with the PHERAstar FS.



**Fig. 5:** HTRF<sup>®</sup> ratios obtained for the IP-One assay with a CCD-based HTS Device.

**Table 1:** Speed and assay quality are compared between the PHERAstar FS and the CCD camera based HTS reader.

	PHERAstar FS	CCD based HTS reader
Read Times (1536)	53 sec	2:12 min
Assay Window	6:1	2:1
Delta F%	490	76
Z' Value	0.70	0.24

## Conclusion

The IP-One HTRF<sup>®</sup> assay from Cisbio was performed on two different HTS microplate readers with different detection technologies. Low affinity compounds which were not discovered with a leading HTS CCD camera based imaging microplate reader were readily resolved with the PMT based PHERAstar FS. This next-generation HTS reader, the PHERAstar FS from BMG LABTECH, represents a new choice for HTS screening assays.



**PHERAstar<sup>®</sup> FSX**

\*The PHERAstar FSX is the newest PHERAstar reader.