

Development of an HTRF-based High Throughput assay for the identification of natural ligands for AMP-activated protein kinase

Yann Ratinaud, Laurent Bultot, Loraine Merminod, Kei Sakamoto, Denis Barron

Introduction

AMP-activated protein kinase (AMPK) is a central cellular energy sensor and regulator (1). To support the research activity on this target the Natural Bioactives and Screening department has the mission to screen natural products that could positively modulate the activity of this enzyme. A radiometric assay was originally developed for our research purposes. Upon activation, AMPK transfers a radiolabeled γ -phosphate (^{32}P) from an ATP to a positively-charged peptide substrate. The phosphorylated substrate binds to a phosphocellulose paper and the enzyme activity can be evaluated by counting the radioactivity(2). Although very sensitive and accurate this methodology is not suitable for high-throughput screening (HTS) campaigns as it normally requires several manual steps and also consumes/produces a substantial amount of radioactivity and waste materials. In contrast, the Homogenous Time Resolved FRET technology (HTRF) is well adapted to HTS and automation: it helps to design simple assays which are homogeneous and usually only requires a few steps that can easily be automatized(3). Two alternative assays from Cisbio based on this technology were evaluated: the HTRF® Transreener® ADP and the HTRF® KinEASE™ STK-S1, and an assay was developed and validated up to a proof of concept screen which was performed on a set of 1000 samples.

Assay Selection

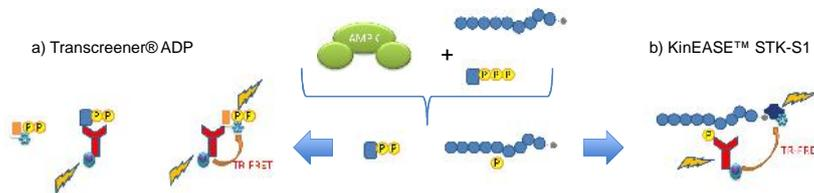


Figure 1: Two distinct approaches based on HTRF technology are used to Monitor AMPK activity: a) Quantification of ADP produced by competition using Transreener® ADP kit; or b) direct quantification of a substrate phosphorylation using KinEASE™ STK-S1 kit

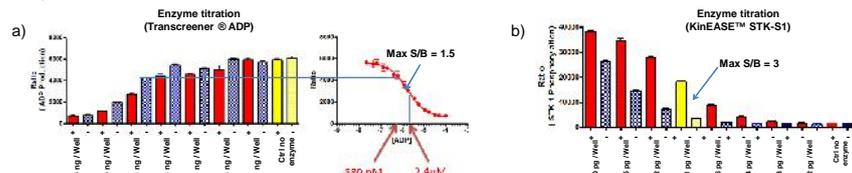


Figure 2: Various concentrations of enzyme are tested in presence (+) or in absence (-) of 30 μM AMP. Reaction is monitored either with a) Transreener® ADP kit, or b) with KinEASE™ STK-S1 kit. The results are reported using the ratio signal (665/620):

Assay development

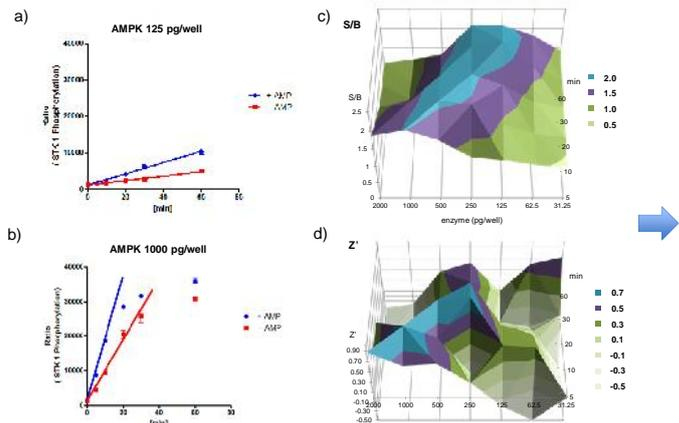


Figure 3: The influence of the reaction duration on substrate phosphorylation is measured in presence or in absence of 30 μM AMP using 125pg/well (a) or 1000 pg/well (b) of AMPK. Signal to background (c) and Z' factor (d) are then calculated on a broader condition test.

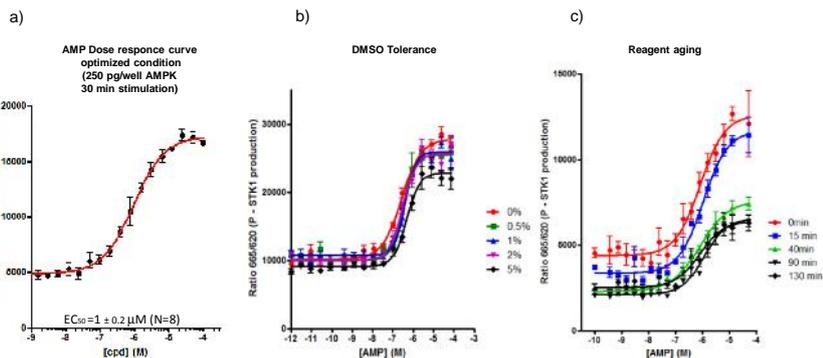


Figure 4: a) AMP titration is performed first in optimized conditions; b) The effect of % DMSO on AMP titration is evaluated; c) Finally the stability at room temperature of the various reagents is estimated by starting the experiment after 5 different waiting times.

Automation



Final Throughput:
- 320 compounds / plates
- 8 plates / run (2 hours)

→ Up to 10 000 samples / day

Figure 5: Automation on the PerkinElmer® Staccato platform. TR-FRET signal is measured off-line using Biotek® SynergyNeo

Proof of concept screening

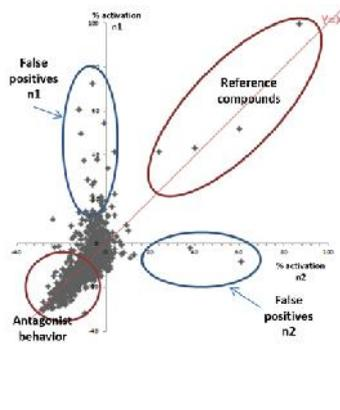


Figure 6: Comparison of n1 Vs n2 % of activation on a proof of concept screening of 1000 samples

Conclusion

- Transreener® ADP kit allows a full flexibility on substrate selection and can be used as a universal readout for all type of kinases
- Direct phosphosubstrate quantification using KinEASE™ STK-S1 requires less enzyme and gives a better S/B ratio (Figure 2)
- The enzyme concentration and the reaction duration are two parameters that greatly influence kinase reaction (Figure 3a,3b)
- Optimal assay conditions can be obtained by co-optimization of enzyme concentration and reaction duration (Figure 3c,3d)
- AMP EC50 determined in the optimized conditions of $1 \pm 0.2 \mu\text{M}$ (Figure 4) matches the one obtained on the original radiometric assay
- DMSO up to 2% has no impact on AMPK enzymatic reaction. Samples will be tested at a final concentration of 1% DMSO.
- After 15 min at room temperature the enzyme activity is reduced. Freshly prepared enzyme must be used for each run on the robotic platform.
- Automation of this assay enabled to reach a throughput of up to 10000 samples/ day (Figure 5).
- Proof of concept screening results from 1000 compounds: (Figure 6)
 - No hits were identified,
 - 14 samples were false positives
 - 17 to 20% of the samples presented an antagonist behavior
 - 5.6 % presented interference with the HTRF technology

References:

1) AMP-activated protein kinase: a target for drugs both ancient and modern. *Chem Biol* 2012 Oct 26;19(10) Hardie DG, Ross FA, Hawley SA.
2) Mechanism of action of compound-13: an 1-selective small molecule activator of AMPK. *Chem Biol* 2014 Jul 17;21(7):866-79. Hunter RW, Forest M, Bultot L, Fullerton MD, Deak M, Ross FA, Hawley SA, Shiro N, Viollet B, Barron D, Kemp BE, Steinberg GR, Hardie DG, Sakamoto K.
3) Development of a HTRF® Kinase Assay for Determination of Syk Activity. *Curr Chem Genomics*. 2008; 1: 20-26. Christopher Harbert, Jeannette Marshall, Sharon Sch, and Krista Steger