



AstraZeneca

Take advantages of HTRF to Simplify HTS Cascade

Jade Zheng, Yue Shao, & Jianwei Liu

Innovation Center China, AstraZeneca, Shanghai, China

Abstract

HTRF assay is homogeneous, very sensitive, and easy to miniaturize. So, it has been used in HTS frequently. Its features of time-resolved and longer emission wavelength, as well as ratio-metric readout result in little potential interference from colored compounds.

Taking these advantages of this assay format, we developed a HTRF assay with high performance for a given target and then simplified the HTS cascade by eliminating single-poke confirmation and the artifact screening.

With this simplified cascade, we delivered final HTS results with excellent outcome in less time and fewer cost.

Platform Validation and HTS

HTS Assay Validation

- ✓ Well2well: ~6% CV
- ✓ Plate2plate & Day2day variations: Reference IC50 < 2x
- ✓ Reference IC50 and S/N (~5): Stable within 24h
- ✓ The 3SD cut-off: ~25%.
- ✓ Double blind test: Captured 40 points of > 30% inhibition with 100% match with the compound map (10 pts of each IC30, 40, 50, 60 in random positions: total 40 points)

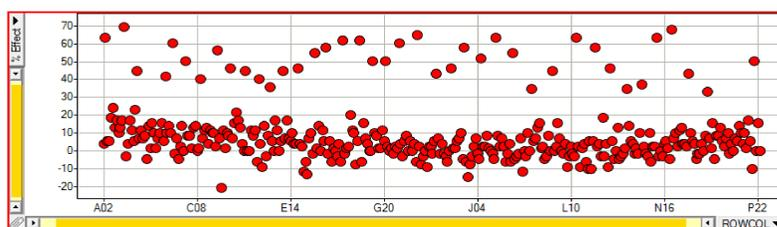


Figure3. Results of double blind test

Normal HTS vs Simplified HTRF HTS

Normal HTS Cascade

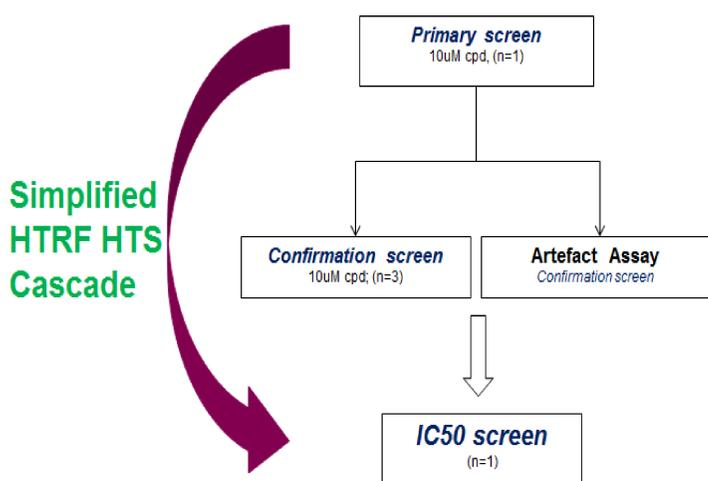


Figure1. Comparison between normal and HTRF HTS cascade

Robust HTS Performance

Validation Run

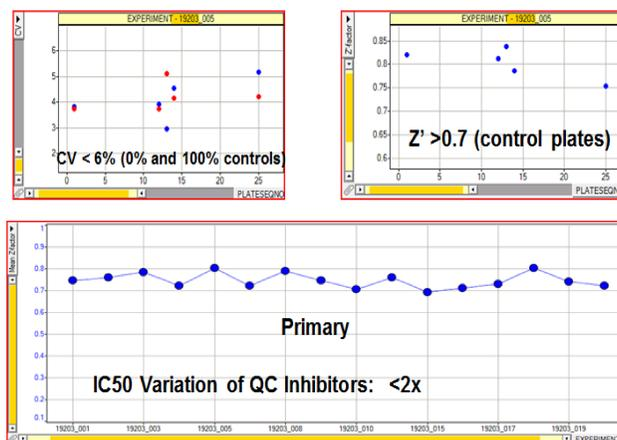


Figure4. Quality Control of HTS performance

Introduction

Assay Description

HTRF KinEASE™ – TK is a generic method for measuring tyrosine kinase activities using one substrate and a universal detection system.

✓ **Enzymatic step:** During this step, the kinase will phosphorylate the substrate. The TK Substrate – biotin is incubated with the kinase. ATP is added to start the enzymatic reaction.

✓ **Detection step:** The detection reagents will catch the phosphorylated substrate. The resulting TR-FRET signal is proportional to the phosphorylation level. The TK-Antibody labeled with Eu3+ - Cryptate and streptavidin – XL665 are then added with EDTA (used to stop the kinase activity).



Step 1: Kinase reaction

- 15 nl compound
- 3 µl Enzyme Mix
- 3 µl TK-substrate

Step 2: Detection

- 6 µl Detection Mix
- Incubation 1h at RT and read on a HTRF reader*

Figure 2. Principle of the HTRF® KinEASE™ - TK assay.

Results of HTS

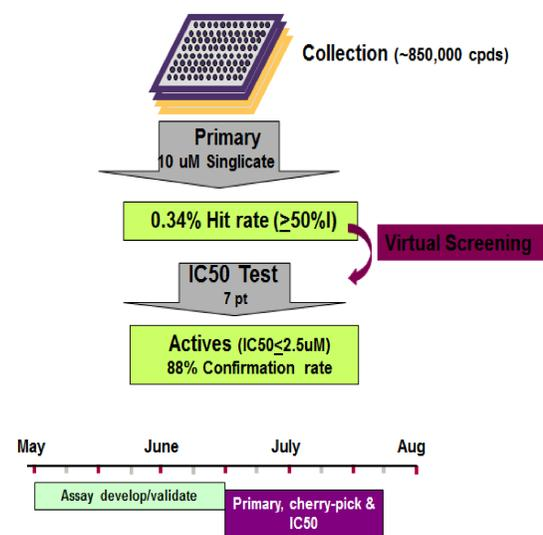


Figure 5. Working model of HTRF HTS

Conclusions

- **HTRF has very low artifact rate:** Artifact screening won't add significant value.
- **HTRF has excellent performance and thus high confirmation rate:** Confirmation screening may not need.
- **Therefore, simple cascade may be applied to HTS using HTRF:** Save time and money.