

Using High Throughput SPA To Screen for CYP2D6 Interactions

A. Hopkins¹, M.J. Price-Jones¹, K.T. Hughes¹, P.J. Eddershaw², S.R. Hood², A.J. Woodrooffe², M.H. Tarbit².

¹ Amersham Biosciences UK Limited, Amersham Place, Little Chalfont, Buckinghamshire, England HP79NA

(Telephone: +44 29 20526417, Fax +44 29 20526474, e.mail Cardiff.SPA.Group@eu.amershambiosciences.com*)

² GlaxoWellcome Research & Development, Technology and Informatics Group, Bioanalysis & Drug Metabolism, Park Road, Ware, England SG12 ODP.

Introduction

Scintillation Proximity Assay (SPA) is an innovative approach to high throughput screening which allows the rapid and sensitive assay of a wide variety of biological processes in a homogeneous system. It is versatile, can be readily automated and assay costs are reduced as the requirement for filters, washing steps and scintillation reagents is eliminated.

The SPA principle is based on the observation that in aqueous solution weakly emitting β -isotopes such as [³H] need to be close to scintillant molecules to produce a light signal. If not, the energy is dissipated into the aqueous solvent.

We have developed a scintillation proximity assay (SPA) using yttrium silicate (YSi) beads, suitable for the measurement of CYP2D6 interactions through the inhibition of [³H]imipramine binding to recombinant CYP2D6 co-expressed with NADPH-cytochrome P450 reductase (CYPOR)⁽¹⁾. Residual binding of the radioactive ligand is determined by scintillation counting following incubation of the test compound with CYP2D6/CYPOR and NADPH (1mM) for up to 30 minutes at 37°C.

P450 SPA competitive binding assay

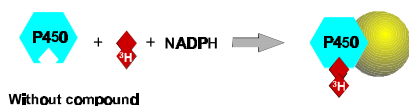


Figure 1. P450 SPA assay concept.

Method

Using a 96-well microplate, 10 μ g CYP2D6 (0.45pmol) in 25 μ l PBS was pre-incubated with the test compound (final concentration 10 μ M) at 37°C for 15 minutes.

This was followed by the addition of 0.1 μ Ci [³H]imipramine (Amersham Biosciences, TRK553, 20-30Ci/mmol, ~150-200,000 liquid scintillation counts added) and 1mM NADPH, 25 μ l each.

The reactions were further incubated for 30 minutes at 37°C. 1mg YSi SPA beads (Amersham Biosciences RPNQ0013) was added to each well, the plate was shaken for 10 minutes and counted using a Wallac MicroBetaTM scintillation counter. Total assay volume was 150 μ l. Results may be obtained for up to 18 hours after the bead addition (see Figure 2).

Results

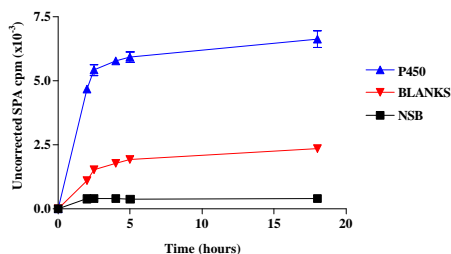


Figure 2. Typical CYP2D6 assay results in the absence of test compounds, shown as means, +/- SEM, (n=3).

"Blanks" are microsomes prepared from mock-infected cells, this signal was subtracted from the P450 counts to give corrected specific SPA cpm. NSB (non-specific binding) was calculated from wells containing PBS, [³H]imipramine and NADPH only.

The uncorrected P450 SPA signal generated was typically >4000cpm. Signal:noise, calculated as the P450 SPA signal divided by NSB, was generally > 10:1.

This assay has been successfully used to distinguish a panel of known compounds that bind to CYP2D6 from a series of P450 probe compounds. The screening results shown as SPA cpm in Table 1 and Figure 3 are from counting the plates immediately, and are means, +/- SEM, (n=3).

| ID | COMPOUND | P450 | BLANKS | CORRT ¹ | % OF CONTROL |
|---------|-----------------|------|--------|--------------------|--------------|
| CONTROL | NO COMPOUND | 5023 | 1100 | 3923 | 100 |
| 1 | BUFURALOL | 1921 | 1139 | 782 | 20 |
| 2 | ONDANSETRON | 2501 | 1152 | 1349 | 34 |
| 3 | AMITRIPTYLINE | 1373 | 1035 | 338 | 9 |
| 4 | DEXTROMETHORPAN | 1365 | 1137 | 228 | 6 |
| 5 | IMIPRAMINE | 1329 | 1098 | 231 | 6 |
| 6 | PROPANALOL | 1520 | 1189 | 331 | 8 |
| 7 | KETOCONAZOLE | 3141 | 1195 | 1946 | 50 |
| 8 | FURAFYLLINE | 4045 | 1322 | 2723 | 69 |
| 9 | PHENACETIN | 5602 | 1252 | 4350 | 111 |
| 10 | CHLORZOXAZONE | 4957 | 1352 | 3605 | 92 |

Table 1 Results and details of the compounds screened at a final concentration of 10 μ M.

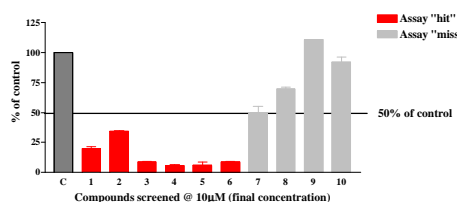


Figure 3 Typical CYP2D6 screening results.

From a panel of 10 P450 probe compounds, 6 were correctly identified as having interactions with CYP2D6.

This assay can be used to generate IC₅₀ values (as shown in Figure 4 below) with, for example, quinidine, a well known inhibitor of CYP2D6.

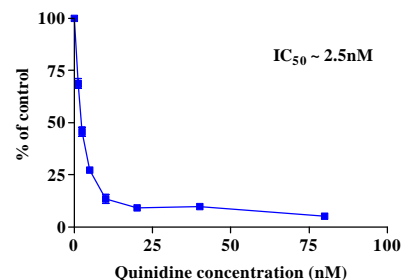


Figure 4 Typical IC₅₀ result achieved with quinidine, a potent inhibitor of CYP2D6.

CONCLUSIONS

- We have developed an SPA assay for the measurement of CYP2D6 interactions which is suitable for the rapid screening of compounds in microplate format.
- The assay utilises [³H]imipramine and has been successfully used to distinguish known compounds that bind to CYP2D6 from compounds that bind to other CYP isoforms.
- An IC₅₀ value of ~2.5nM was generated with quinidine, a well known CYP2D6 inhibitor.

Reference

- HOOD, S.R. *et al.*, *Methods in Mol. Biol.*, **107**, 203-218, (1998).

This poster has been produced as a result of a Technology Transfer collaboration between: