

# A HIGH THROUGHPUT SCREENING ASSAY FOR THIAMINE PYROPHOSPHATE KINASE ACTIVITY USING SPA

<sup>1</sup>Greet Vanhoof, <sup>2\*</sup>Jenny Berry, <sup>2</sup>Mark Harvey, <sup>2</sup>Molly Price-Jones and <sup>2</sup>Kelvin Hughes.

<sup>1</sup>Janssen Research Foundation, Beerse, Belgium.

<sup>2</sup>Amersham Biosciences UK Limited, Amersham Place, Little Chalfont, Buckinghamshire, England, HP7 9NA (Telephone +44 (0) 29 2052 6417, Fax +44 (0) 29 2052 6474, e: mail Cardiff.SPAGroup@eu.apbiotech.com)

## Introduction

Thiamine pyrophosphate kinase (TPPK: E.C. 2.7.6.2) converts thiamine to thiamine pyrophosphate (TPP) which acts as the major metabolic thiamine signal within the cell<sup>(1)</sup>.

TPP functions as a cofactor for many enzymes involved in carbohydrate metabolism, and in *Schizosaccharomyces pombe*, has been shown to be a critical metabolic signal for mating. TPP is also thought to be important in phosphate metabolism and growth in yeast<sup>(2)</sup>. Therefore, inhibition of TPPK may be a target for anti-fungal therapy.

Current methods for analysing TPPK activity such as HPLC<sup>(3)</sup> and ion exchange chromatography<sup>(4)</sup> are not amenable to high throughput screening; therefore, we have developed a scintillation proximity assay (SPA) to measure TPPK activity.

In this assay, TPPK from *Saccharomyces cerevisiae* was used to convert [<sup>3</sup>H]thiamine to [<sup>3</sup>H]TPP which was selectively captured using yttrium silicate (YSi) SPA beads. (Fig. 1).

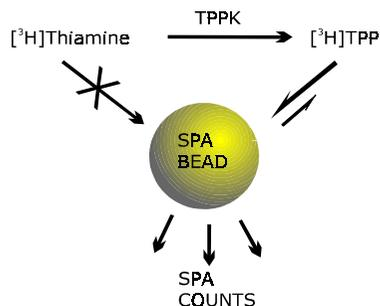


Figure 1. Diagrammatic representation of the TPPK SPA assay.

## Method and Results

SPA assays contained TPPK (0.5µg), unlabelled thiamine (10nmol) and [<sup>3</sup>H]thiamine (70pmol, 1µCi) in a total volume of 100µl 5mM glycine buffer, pH 8.6, containing 13µM ATP and 2mM MnSO<sub>4</sub>. Assays were incubated at room temperature for 20 minutes before being stopped by the addition of 50µl de-ionized water containing 1mg YSi SPA beads. Non-specific binding was determined in the absence of TPPK.

Assays were counted for 1 minute/well using a TopCount™ microplate scintillation counter. Library compounds were added in DMSO to give a final concentration of 1% DMSO (v/v) and 10<sup>-5</sup>M test compound in the assay well.

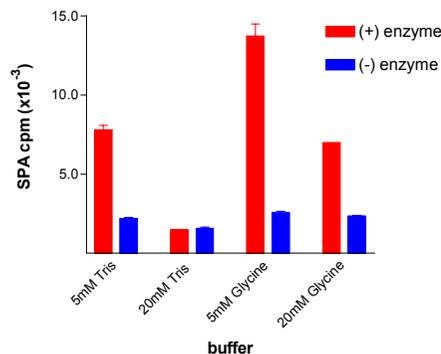


Figure 2. Effect of buffer type and concentration on SPA assay signal. Both tris and glycine buffers (containing 13µM ATP and 2mM MnSO<sub>4</sub>) were used at pH 8.6. Values are means ±SEM (n=2).

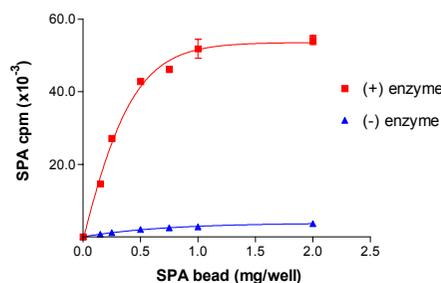


Figure 3. Effect of increasing concentrations of YSi bead on SPA assay signal. Assays contained 1.0µg enzyme/well. Values are means ±SEM (n=2).

As can be seen from Figure 2, both the type of buffer, and the buffer concentration, appeared to have a considerable effect on the SPA cpm obtained. The use of glycine buffer resulted in ~40% increase in SPA cpm compared to tris buffer. Increasing the glycine buffer concentration from 5 to 20mM appeared to reduce the SPA cpm by ~50%.

A titration of SPA bead (Fig.3) indicated that the maximum signal: background was obtained with 1mg YSi bead/well.

Experiments indicated that the assay signal was linear up to 1.25µg/well added enzyme (Fig.4).

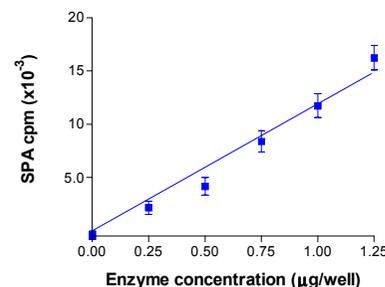


Figure 4. Effect of increasing enzyme concentration. Values are means ±SEM (n=3).

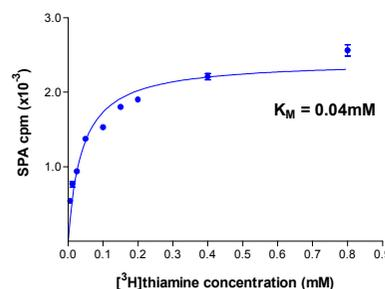


Figure 5. Michaelis-Menten plot for TPPK with [<sup>3</sup>H]thiamine as substrate. Values are means ±SEM (n=3).

Non-linear regression analysis of the substrate concentration curve obtained with [<sup>3</sup>H]thiamine (Fig.5) gave a K<sub>M</sub> value of 40µM (± 8µM). This is in good agreement with the previously reported K<sub>M</sub> values for prokaryotic TPPK (75µM for *Saccharomyces cerevisiae* TPPK<sup>(5)</sup>, 6µM for *Saccharomyces pombe* TPPK<sup>(2)</sup> and 38µM for *Paracoccus denitrificans* TPPK<sup>(6)</sup>).

Pyriothiamine, a substrate analogue and thiamine antagonist, is thought to be a direct competitive inhibitor of TPPK. Reported K<sub>i</sub> values for pyriothiamine of the prokaryotic TPPK are 6µM<sup>(2)</sup> and 19µM<sup>(6)</sup>. As can be seen from Fig.6, inhibition studies indicate that pyriothiamine does inhibit *Saccharomyces cerevisiae* TPPK with an estimated IC<sub>50</sub> value of 19µM.

The robustness of this SPA assay was tested using a high throughput screening system, screening 42 x 96-well plates of our compound library, where each plate

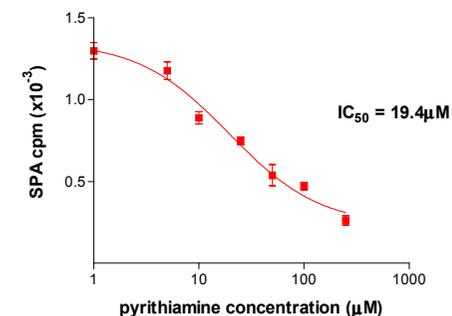


Figure 6. Inhibition of TPPK by pyriothiamine. Values are means ±SEM (n=3).

included 8 wells of positive controls (in the presence of TPPK enzyme) and 8 wells of negative controls (in the absence of TPPK), testing a total of 3360 compounds. The average of the cpm values for the positive controls was 6245 cpm with a CV of 7%. The average cpm value for the negative control was 816 cpm with a CV of 8%. To evaluate the quality of the HTS we calculated the Z' value<sup>(7)</sup> and found it was 0.7. The quality of the assay allows implementation in an HTS system with delivery of precise and reproducible data.

## Conclusions

- We have developed a high throughput screening assay to measure the inhibition of TPPK activity.
- The assay is robust, and can be used for automated screening of a large number of compounds for discovery of hits and development of therapeutic leads.

## References

1. VOSKOBOYEV, A.I., et al., *Ann. N. Y. Acad. Sci.*, **40**, 161-76 (1982)
2. FANKHAUSER, H., et al., *J. Biol. Chem.*, **270** (47) 28457-62 (1995)
3. EGORAMAIPHOL, S., et al., *Int. J. Vitam. Nutr. Res.*, **61** (4) 334-8 (1991)
4. DEUS, B., et al., *Methods Enzymol.*, **62**, 103-4, (1979)
5. KAZIRO, J. et al., *J. Biochem.*, **46**,1523-39 (1959)
6. SANEMORI, H. et al., *J. Biochem.*, **88**, 223-30 (1980)
7. ZHANG, et al., *J. Biomolec. Screening*, **4**, 67-73 (1999)