

THYMIDINE KINASE 1 IN THE STUDY OF SOLID TUMORS

Thymidine Kinase 1 (TK1) has been long known as a valuable biomarker of cellular proliferation. However, previous methods have been based on enzyme activity measurements that may be subject to interference and may underestimate the forms of TK1 found in the serum of subjects with solid tumors.

The AroCell TK 210 ELISA kit is a new valuable biomarker for studying the rate of proliferation and cell turnover of solid tumors.

It brings the specificity of immunoassay to this field and offers improved accuracy when studying serum TK1 derived from patients with solid tumors.

The AroCell TK 210 ELISA kit will provide new opportunities for studying cellular proliferation, tumor cell turnover and therapy response in subjects with solid tumors.

Thymidine Kinase 1 in Studying Solid Tumors

Tissue TK1

DNA replication requires the production of phosphorylated thymidine. There are two enzymes that can perform this, Thymidine Kinase 2 (TK2), found in mitochondria, which is constitutively expressed and Thymidine Kinase 1 (TK1) the expression of which varies during the cell cycle. TK1 concentrations in the cell are low in the G0/G1 phase (resting phase) of the cell cycle but increase during the S/G2 phases when DNA synthesis occurs and then decrease during mitosis¹. The presence of TK1 in cells is an indicator of active cellular proliferation.

TK1 up-regulation may be an early event in cancer development and TK1 may even be elevated in pre-cancerous conditions². Increased TK1 expression is often associated with increased expression of cell proliferation markers such as the Ki-67 antigen and proliferating cell nuclear antigen (PCNA)^{3,4,5} although studies have shown that TK1 may be a more useful as a proliferation marker than either of them⁶. TK1 has the further practical advantage in that it is measurable in both tissues and serum, simplifying serial testing.

Increased cellular proliferation is a hallmark of malignancies and TK1 has been found to be significantly over-expressed compared to normal tissue in many solid tumors including lung, colon, prostate, esophagus, stomach, liver, and renal cancers⁷.

Tissue TK1 activity may aid in making prognoses and in the selection of therapeutic alternatives in that subjects with high tissue levels may have a shorter survival time and that subjects with lower levels may respond better to anti-hormonal, therapy⁸.

TK1 activity in serum

95% of serum TK1 activity in cancer patients seems to be tumor derived, making TK1 an excellent indicator of cellular proliferation⁹ and a useful complement to immunohistological testing for proliferation biomarkers. TK1 enzyme activity has been shown to be elevated in subjects with many forms of cancer, including leukaemia, lymphoma, prostate, breast, lung, sarcoma and colon cancer patients¹. Elevations in serum TK1 may be an early event and subjects can have elevated serum TK1 at presentation and high pre-treatment TK1 activities have been associated with shorter progression-free and overall survival in subjects with breast cancer¹⁰. Conversely, subjects with cancer but lower serum TK1 values may have improved survival¹¹. Continued elevations in serum TK following surgery may indicate the residual tumor and an increased likelihood of relapse¹².

TK1 serum activity is a valuable biomarker but, being an enzyme activity assay, it has limitations regarding specificity and sensitivity in that serum contains inhibitors of TK1 activity and that serum contains multiple forms of TK1 with differing specific activities.

The development of immunoassays for TK1 protein mass can reduce these problems and increase the diagnostic value of TK1 as a biomarker.

TK 210 as a cell proliferation biomarker

As mentioned above, although TK1 activity is a valuable biomarker it has limitations in that most current methods are based on enzyme activity assays¹³. The AroCell TK 210 ELISA kit, based on two monoclonal antibodies specific for a unique TK1 amino acid sequence (194-225), brings the specificity, sensitivity and robustness of immunoassay to the assay of serum TK1¹⁴.

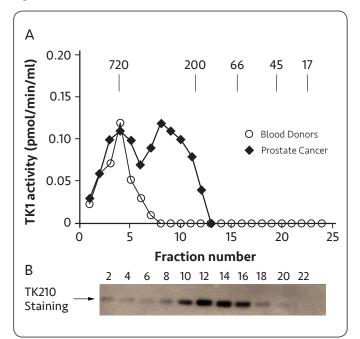
Studies at AroCell demonstrated that certain TK1 forms that can be found in subjects with solid tumors differ from those found in normal serum and hematological malignancies. These forms differed in both molecular weight and specific activity and by also measuring these forms the TK 210 immunoassay offered greater discrimination and sensitivity. Furthermore, the use of ELISA simplifies the assay procedure and avoids the effects of interfering substances.

The AroCell TK 210 ELISA kit enables TK1 protein levels in subjects with solid tumors to be more clearly distinguished from those found in healthy subjects.

Analysis of Serum TK1 Protein Levels with TK 210-Immunoassay

Greater sensitivity and specificity for tumor associated TK1

Figure 1



Tissue TK1 is formed of four identical 25kDa units but serum TK1 enzyme activity is found associated with proteins covering a wide range of molecular weights and this distribution differs between healthy subjects and those with malignancies¹⁵. Figure 1A shows how the distribution of TK1 activity in serum from a subject with prostate cancer differs from that in serum from a healthy blood donor. In the blood donor serum, TK1 activity was found as a single peak of MW between 500-720 kDa whereas in prostate cancer sera, TK1 activity eluted as a broad peak in fractions corresponding to the MW range 200-720 kDa.

Western blot analysis of serum from a prostate cancer subject (figure 1B), using the TK 210 antibody as a probe, showed that TK1 units (25 kDa) were found in fractions corresponding to a wide range of molecular weights, especially below 500 kDa where enzyme activity was low or absent from blood donor sera. Furthermore, there are low molecular weight TK1 fractions (< 200kDa) in serum from prostate subjects where TK1 activity is totally lacking (fractions 14 onward in figure 1B). These results, and similar data from breast cancer, indicate that there are TK1 protein fractions in serum from subjects with solid malignancies which are enzymatically inactive but can be detected and measured immunologically.

TK 210 is more discriminating in solid tumors

TK 210 concentration and TK1 activity in breast cancer

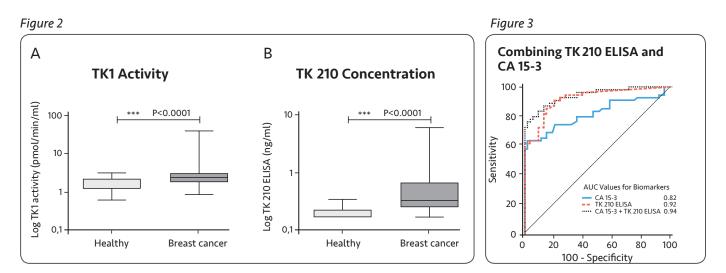


Figure 2 compares TK1 activity levels and TK 210 levels in subjects with breast cancer and blood donors. Note that mean TK 210 ELISA levels differ between breast cancer subjects and blood donors by a factor of five, while TK1 activity levels differ by only a factor of three. Comparing the TK1 activity and TK 210 ELISA as Receiver Operator Characteristic (ROC) curves (Figure 3) shows the greater diagnostic accuracy of TK 210 ELISA (Area under the curve (AUC)

0.92 v 0.82). Interestingly, both showed greater power than the classic tumor biomarker CA15-3 (0.82). Combining TK 210 ELISA and CA15-3 increased the diagnostic power even further to give an AUC value of 0.94. Sensitivity was increased to 0.76 while specificity was retained (0.98). This shows the potential additive value of TK 210 to existing biomarker panels¹⁴.

AroCell TK 210 ELISA as a biomarker for solid tumors

- TK1 is an early biomarker of proliferation
- Serum levels reflect tumor proliferation
- TK1 can be used to study many tumors
- TK1 may aid in prognoses and therapy selection
- Better discrimination between healthy controls and subjects with solid tumors
- Unaffected by serum inhibitors
- Translational biomarker
- Standard ELISA procedure
- CE marked

References

- Topolcan, O. and Holubec, L. (2008). The Role of Thymidine Kinase in Cancer Diseases. Expert Opinion Medical Diagnosis 2(2): 129-141.
- Alegre, M.M. *et al.* (2012). Thymidine Kinase 1 Upregulation Is an Early Event in Breast Tumor Formation. Journal of Oncology. Volume 2012, 5 pages.
- Gasparri, F. *et al.* (2009).Thymidine kinase 1 expression defines an activated G1 state of the cell cycle as revealed with site-specific antibodies and ArrayScan assays. Eur J Cell Biol. 88(12): 779-85.
- Wu, J. *et al.* (2000). A new cell proliferating marker: cytosolic thymidine kinase as compared to proliferating cell nuclear antigen in patients with colorectal carcinoma. Anticancer Res. 20(6C): 4815-20.
- Brockenbrough, J.S. et al. (2009). Thymidine Kinase 1 and Thymidine Phosphorylase Expression in Non–Small-cell Lung Carcinoma in Relation to Angiogenesis and Proliferation. Journal of Histochemistry & Cytochemistry. 57(11): 1087–1097.
- Foekens, J.A. *et al.* (2001). Thymidine Kinase and Thymidylate Synthase in Advanced Breast Cancer: Response to Tamoxifen and Chemotherapy. Cancer Res; 61 (4): 1421-5.
- Alegre, M.M. et al. (2013). Thymidine Kinase 1: A Universal Marker for Cancer. Cancer and Clinical Oncology. 2(1): 2013.
- Guan, H. *et al.* (2009). Thymidine kinase 1 expression in atypical ductal hyperplasia significantly differs from usual ductal hyperplasia and ductal carcinoma in situ: A useful tool in tumor therapy management. Mol Med Rep. 2(6): 923-9.

- Wu, C. *et al.* (2003). Production and characterization of a novel chicken IgY antibody raised against C-terminal peptide from human thymidine kinase 1. J Immunol Methods **277**: 157-169.
- Bjöhle J. et al. (2013). Serum thymidine kinase activity compared with CA 15-3 in locally advanced and metastatic breast cancer within a randomized trial. Breast Cancer Res Treat. 139(3): 751-8.
- Chen, F. *et al.* (2013). Serum thymidine kinase 1 levels predict cancer free survival following neoadjuvant, surgical and adjuvant treatment of patients with locally advanced breast cancer. Molecular and Clinical Oncology. 1: 894-902.
- He, Q. *et al.* (2006). Thymidine Kinase 1 in Serum Predicts Increased Risk of Distant or Loco-regional Recurrence Following Surgery in Patients with Early Breast Cancer. Anticancer Research 26: 4753-4760.
- He, Q *et al.* (2005), Concentration of Thymidine Kinase 1 in Serum (S-TK1) is a More Sensitive Proliferation Marker in Human Solid Tumors Than its Activity. Oncol Rep. **14(4)**: 1013-9.
- Kumar, J.K. *et al.* (2016). A Clinical Evaluation of the TK 210 ELISA in Sera from Breast Cancer Patients Demonstrates High Sensitivity and Specificity at All Stages of the Disease. Tumor Biology. **37**: 11937-11945.
- Jagarlamudi. K.K. et al. (2015). Breast and Prostate Cancer Patients Differ Significantly in Their Serum Thymidine Kinase 1 (TK1) Specific Activities Compared with Those Found in Hematological Malignancies and Blood Donors: Implications of Using Serum TK1 as a Biomarker. BMC Cancer 15:66.

AroCell TK 210 ELISA is for Research Use Only in the USA. Not for use in diagnostic procedures



©2017 AroCell AB. All rights reserved. 16-1MA004B1-04

AroCell AB (publ)., Uppsala Business Park, Virdings allé 32 B, SE-754 50 Uppsala, Sweden e-mail: info@arocell.com, www.arocell.com