

Development of a pan-cancer NGS assay for detection of tumor mutational burden and targeted biomarkers from FFPE samples

Dinesh Cyanam, Vinay Mittal, Nickolay Khazanov, Paul Williams, Santhoshi Bandla, Janice Au-Young, Gary Bee, Sameh El- Difrawy, Aren Ewing, Jennifer Kilzer, Anelia Kraltcheva, Scott Myrand, Yu-Ting Tseng, Warren Tom, Cristina Van Loy, Elaine Wong-Ho, Chenchen Yang and Seth Sadis. Clinical Sequencing Division, Thermo Fisher Scientific, Ann Arbor, MI; Carlsbad, CA; South San Francisco, CA

INTRODUCTION

Next-generation sequencing (NGS) is used to support routine clinical research in oncology with a primary focus on evaluating known oncogenic variants. Effective solutions such as targeted NGS assays allow assessment of hundreds of cancer-related genes simultaneously. Although, the primary focus of targeted assays has been to evaluate known oncogenic variants, the advent of cancer immunotherapies requires that clinical research solutions must also address biomarkers such as Tumor Mutational Burden (TMB) and Microsatellite Instability (MSI) for immune checkpoint inhibitors.

In recent years, TMB has emerged as important biomarker for immunotherapy¹. Several published studies have demonstrated that high TMB is associated with positive response to various immune checkpoint inhibitors.

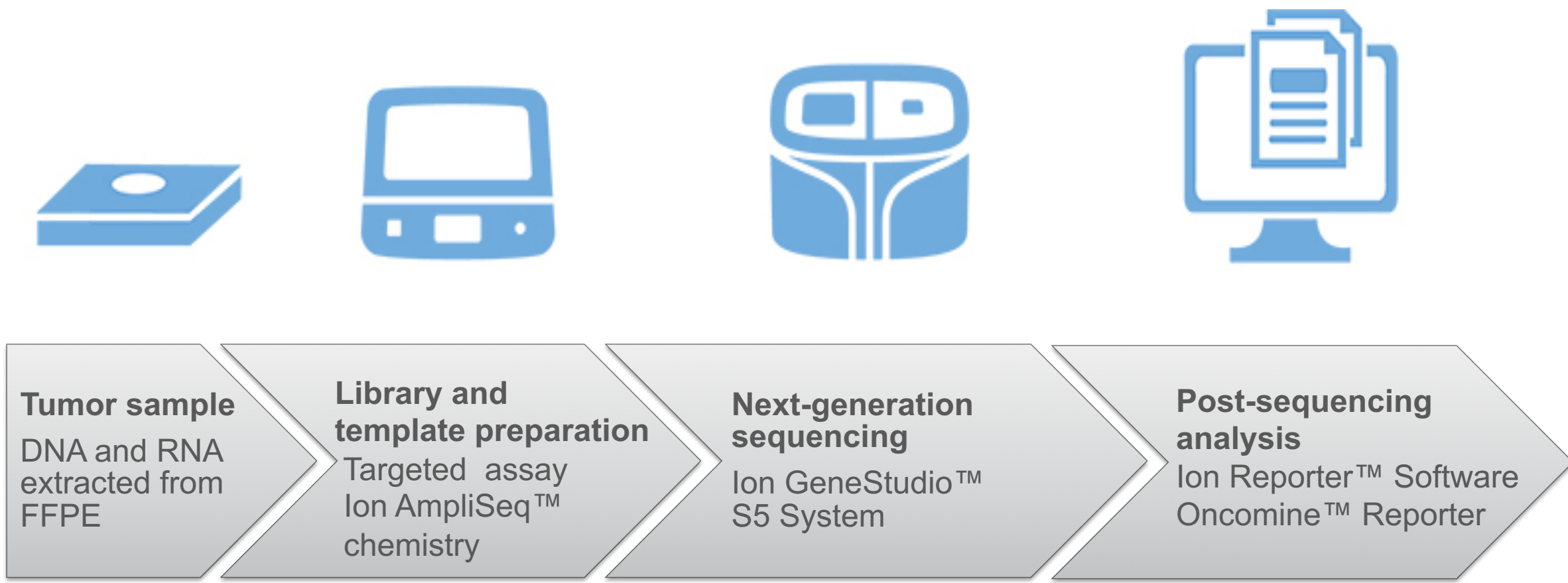
Although TMB can be accurately computed using whole exome sequencing (WES), associated challenges such as high cost, requirement for high starting FFPE material, requirement of tumor matched normal or control samples and limited sample to report software solutions has hindered the adoption of TMB biomarker testing in routine clinical research.

In order to overcome these challenges, we have developed a unified, yet simple multiplex PCR-based target enrichment NGS assay. The assay covers comprehensive targets that are relevant in cancer, has sensitive and specific chemistry to maximize low quantity FFPE tissues, and an automated sample-to-report workflow, that holistically provides an assessment of important cancer biomarkers, including TMB, in a time sensitive manner.

METHODS

Gene content was prioritized based on the relevance and variant prevalence of biomarkers in solid tumors. Additional genomic regions were added to supplement the coding sequence footprint to support TMB. The assay used Ion AmpliSeq™ technology with automated templating on the Ion Chef™ system and sequencing on the Ion GeneStudio™ S5 sequencing platform. An automated tumor-only workflow for variant calling, TMB and MSI estimation and sample quality reporting was provided within Ion Reporter Software. Streamlined access to decision support software is enabled by OncoPrint™ Reporter².

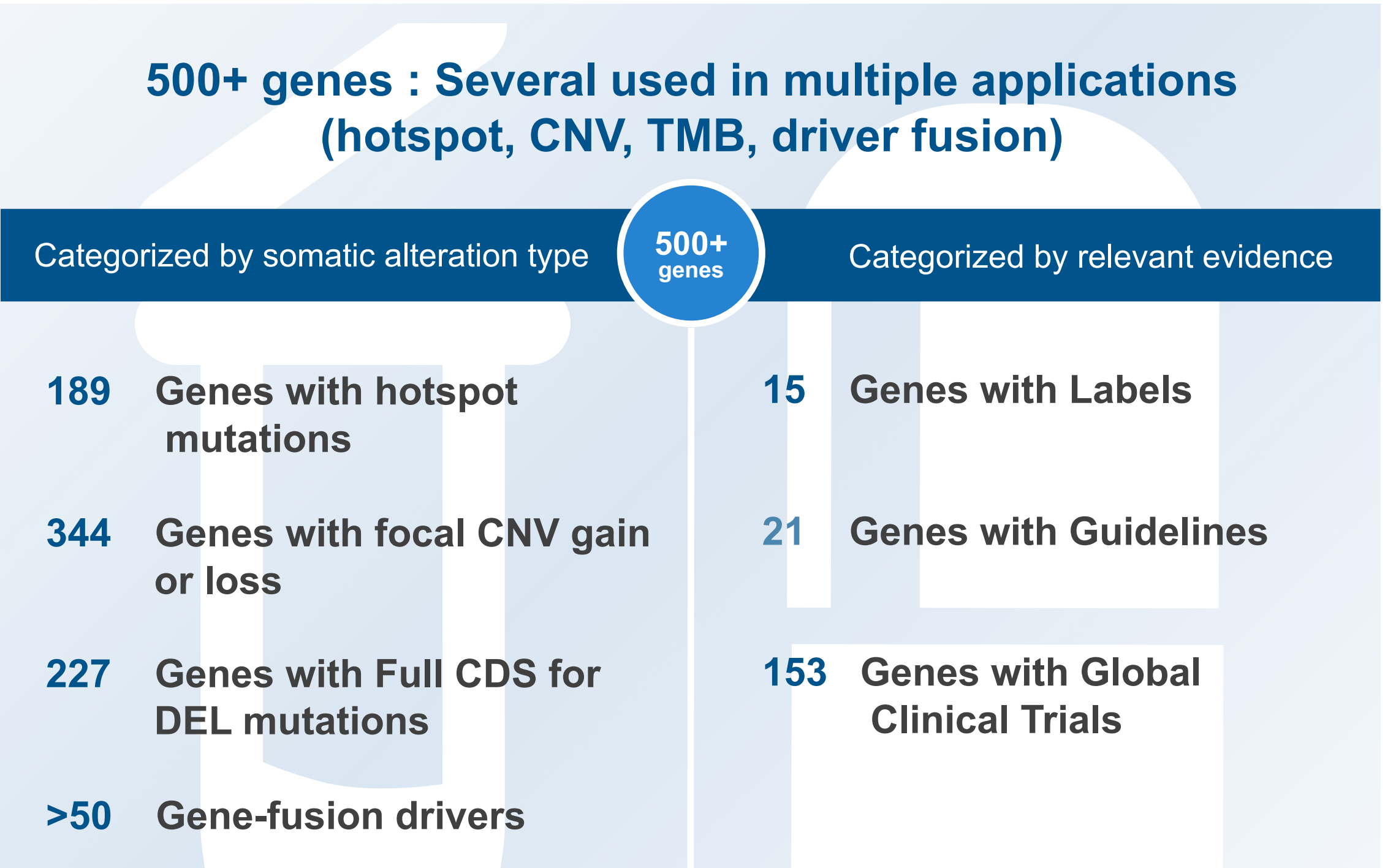
Figure 1. Schematic flow-diagram of the complete workflow



RESULTS

Over 500 genes with DNA based alterations and over 50 RNA fusion drivers are included. More than 13,000 DNA amplicons cover a comprehensive genomic targeted region with a large (>1 MB) coding sequence (CDS) footprint to support high-confidence TMB. FFPE tumor samples from a variety of tissue types were sequenced using the assay. The assay displays high (>95%) uniformity and consistent read depth (>2200x) to support robust variant calling at low allele frequency. In-silico assessment of TMB using publicly available whole-exome cancer sequencing data resulted in high correlation ($R^2 > 0.90$, 0-40 mut/mb) in pan-cancer and specific cancer types including lung, colorectal and melanoma.

Figure 2. Summary of the assay content

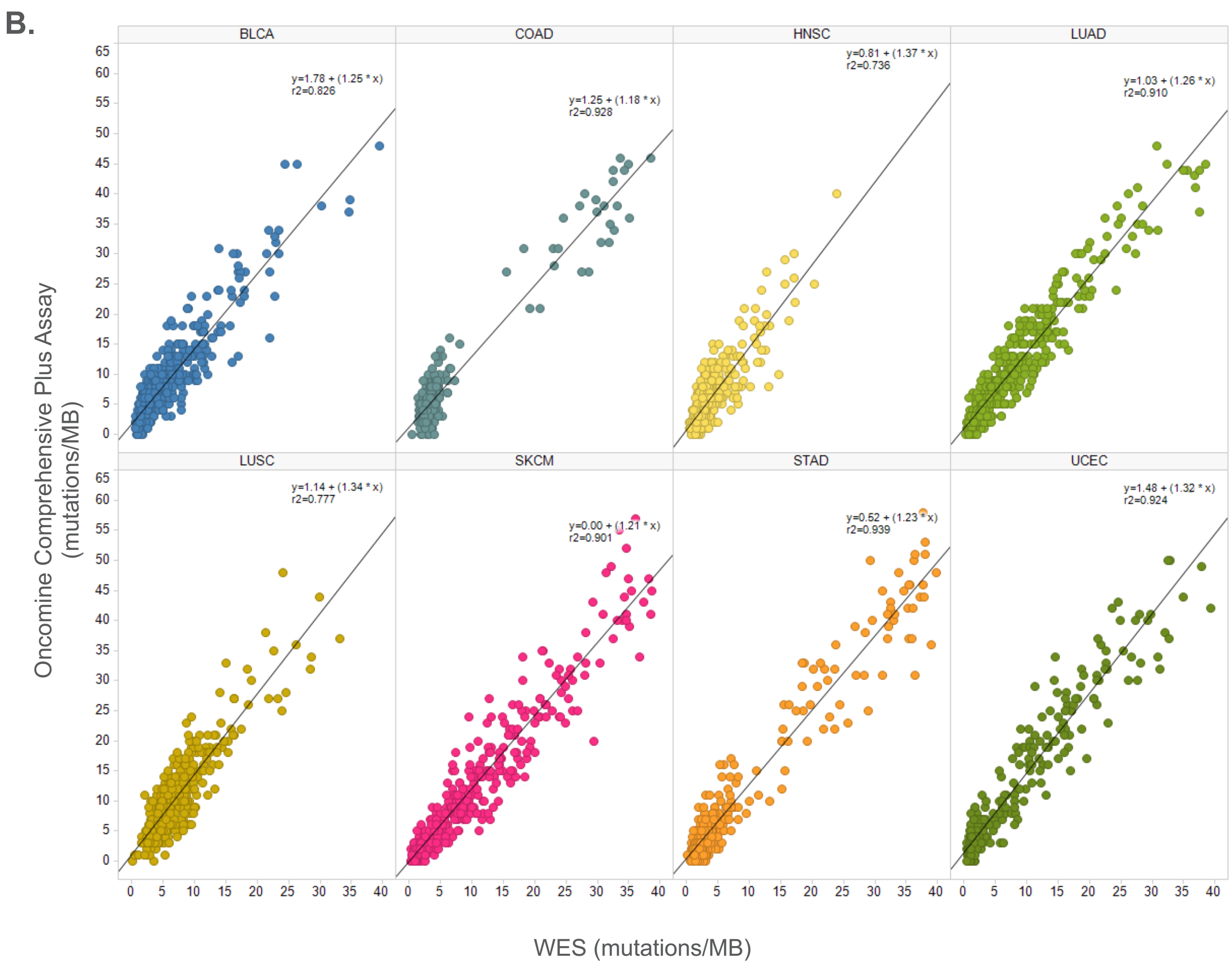
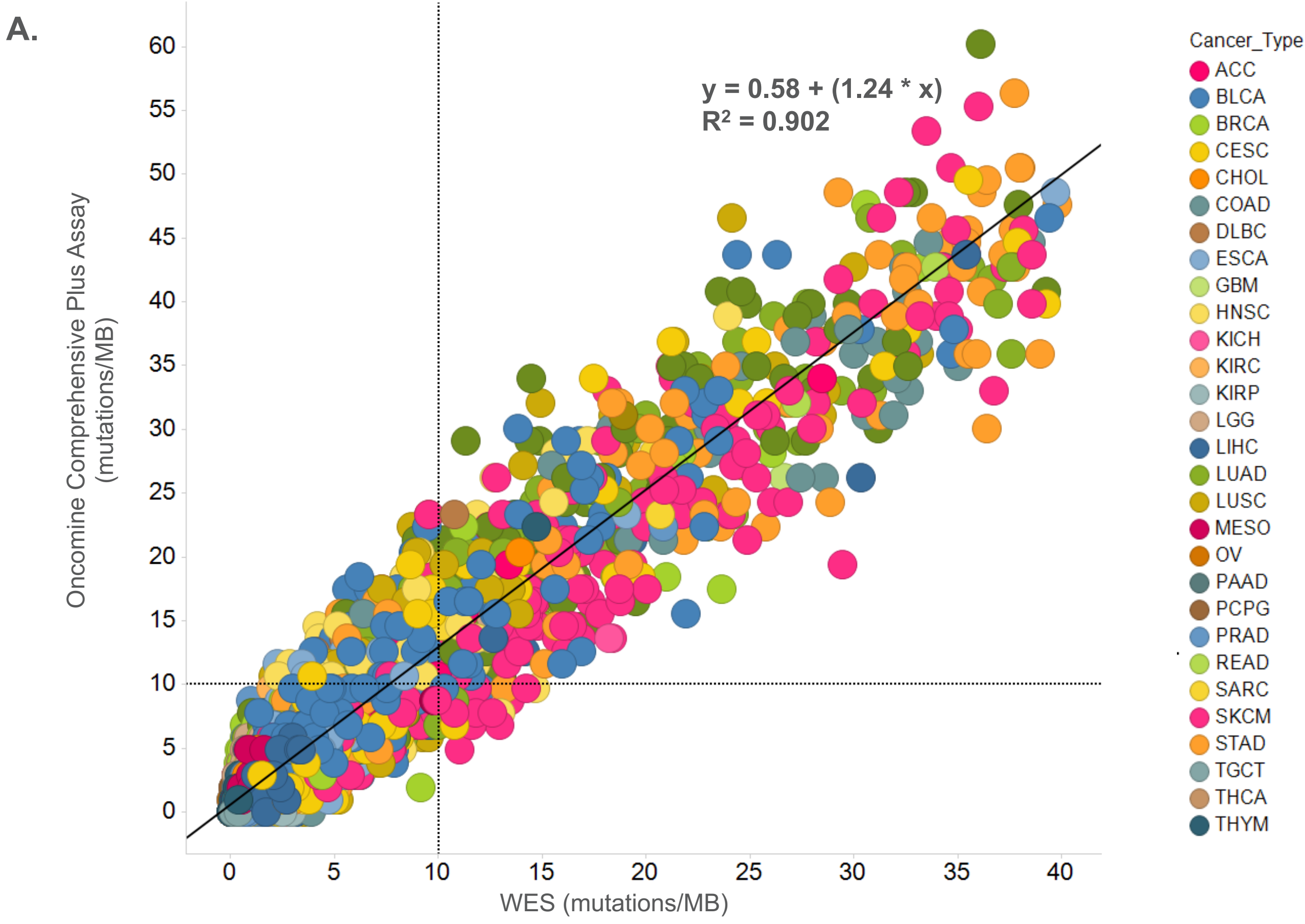


Out of the DNA genes, 189 genes cover important cancer hotspots, 344 cover copy number variants (CNV), while 227 genes have full coding sequence (CDS) coverage for detection of deleterious variants. In addition to fusion drivers, this assay also includes seven genes with intragenic variants. The assay was also designed to maximize the genomic footprint to support high-confidence TMB estimation. The total genomic coverage of the assay is 1.50MB with 1.03MB of coding sequence.

Table 1. OncoPrint Comprehensive Plus Assay Cancer Gene Targets

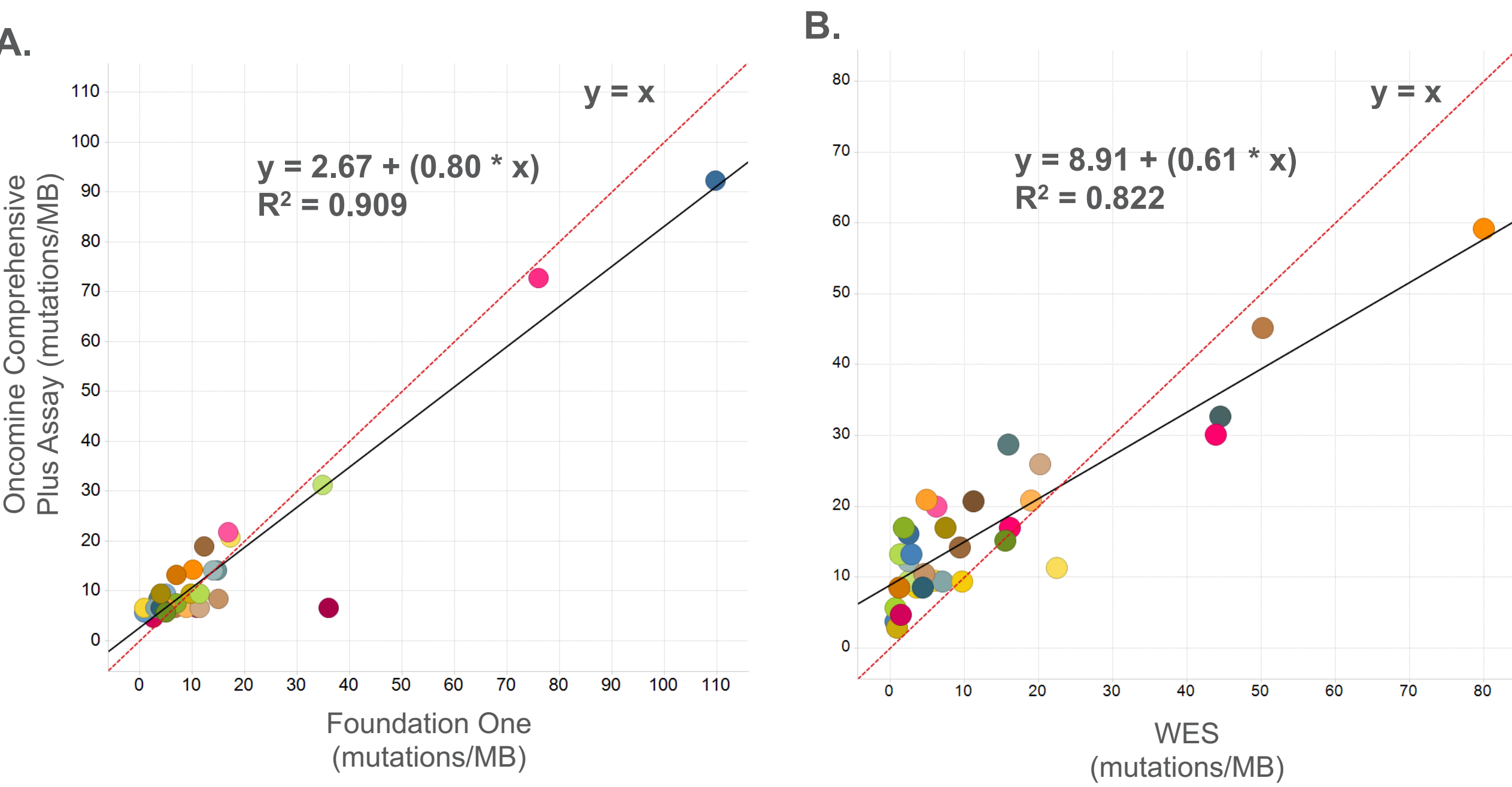
[illegible]

Figure 3. In-silico comparison of the OncoPrint Comprehensive Plus Assay with Whole Exome Sequencing (WES)



Scatter plots (Fig. 3) showing the correlation between the targeted assay (y-axis) and WES (x-axis) mutation counts. WES data was downloaded from TCGA MC33. In-silico analysis was performed to characterize TMB performance of the targeted sequencing assay. Rate of nonsynonymous somatic mutations was computed for WES TMB. Mutations were limited to the targeted assay for predicted TMB. WES TMB strongly correlated ($R^2 > 0.9$) with the assay TMB in pan-cancer analysis (Fig. 3A). Figure 3B displays similar distribution for selected cancer types. Strong correlation ($R^2 > 0.9$) were observed for various cancer types with variations in Bladder Urothelial Carcinoma (BLCA) and Lung squamous cell carcinoma (LUSC).

Figure 4. Comparison of the OncoPrint Comprehensive Plus Assay with Foundation One and Whole Exome Sequencing



Foundation plots (Fig. 4) showing the correlation between the OncoPrint Comprehensive Plus Assay and the Foundation One (Fig 4A) and Whole Exome Sequencing (Fig 4B). OCA Plus Assay shows strong correlation ($R^2 > 0.9$) with Foundation One panel (Fig 4A). For comparison with WES, 35 FFPE tumor samples with matched normal were sequenced on the Illumina platform and the TMB scores were determined. The matched tumor samples were then sequenced on the OncoPrint Comprehensive Plus Assay using IonTorrent sequencing technology and TMB scored were determined. These scores were then compared with the WES TMB scores (Fig 4B).

CONCLUSIONS

A targeted NGS assay was developed to support comprehensive genomic profiling and routine clinical research in oncology that includes endpoints for TMB and MSI. The design and informatics workflow support characterization of mutational signatures and provides normalized TMB estimates with a strong correlation with orthogonal methods. The assay also supports detecting relevant RNA structural alterations from solid tumor FFPEs. Minimal input material requirement and rapid sample to report time will have a high impact on clinical research.

REFERENCES

1. Hellmann et al. Nolumab plus Ipilimumab in Lung Cancer with a High Tumor Mutational Burden. *NEJM* 2018
2. OncoPrime Reporter:
<https://www.thermofisher.com/order/catalog/product/A34298>
3. Ellrott et al. Scalable Open Science Approach for Mutation Calling of Tumor Exomes Using Multiple Genomic Pipelines. *Cell Systems* 2018