

Performance of a Combined SNP and Copy Number Pharmacogenomic Research Panel to Determine Phenotypes of Drug Metabolism Genes

Tibor Fule¹, Kathleen Hayashibara², Pius Brzoska²

¹Thermo Fisher Scientific – Budapest, Hungary, ²Thermo Fisher Scientific – South San Francisco, CA, USA.

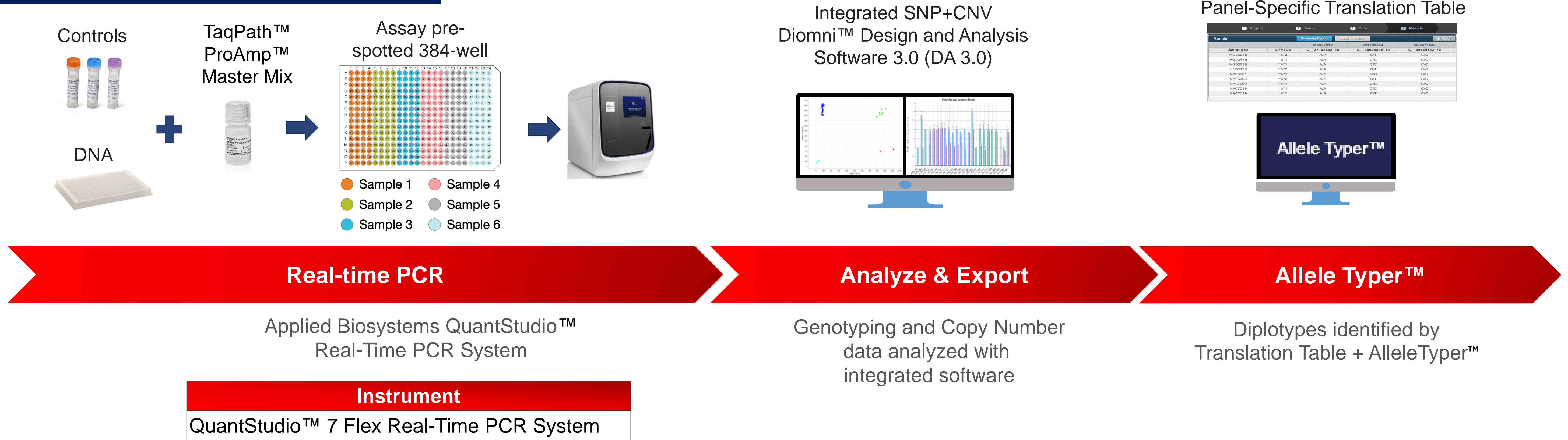
Introduction

Understanding sequence and copy number variations in genes coding for metabolizing enzymes are important when studying drug metabolism. We created a 384-well qPCR research panel containing both TaqMan® SNP Genotyping and Copy Number assays targeting CYP2B6, CYP2C9, CYP2C19, CYP2D6, CYP3A4, CYP3A5, CYP4F2, DPYD, NUDT15, SLCO1B1, TPMT, and VKORC1. Accuracy of phenotypes is presented using Genetic Testing Reference Material (GeT-RM) DNAs containing targeted variants.

Methods

DNAs were normalized to 4 ng/ul, mixed with TaqPath™ ProAmp™ Master Mix, loaded onto pre-spotted assay plates, and run on a QuantStudio™ 7 Flex Real-Time PCR System. CNV and genotyping results were determined by Diomni™ Design and Analysis v3.0 software. Diplotype calls were made using AlleleTyper™ Software v1.0 with translations based on information from PharmGKB.

Product Concept and Workflow



Results

CYP2D6 5'UTR and Exon 9 assays produced 100% call rate and 100% of calls matched consensus results from several platforms. The 56 panel genotyping assays produced results consistent with consensus star alleles in > 95% of sample-target combinations. Sample-target combinations with single diplotype calls matched consensus phenotypes in 100% of cases. Samples with multiple diploypes produced at least one combination matching consensus phenotypes in all cases.

TaqMan SNP genotyping and Copy Number Variation Data

