

A highly specific and sensitive Applied Biosystems™ TaqMan SCID/SMA multiplex assay optimized for dried blood spots

Sonu Baral¹, Kristina Gill¹, Amy Huang¹, Kimitoshi Nakamura², Takaaki Sawada²

BACKGROUND

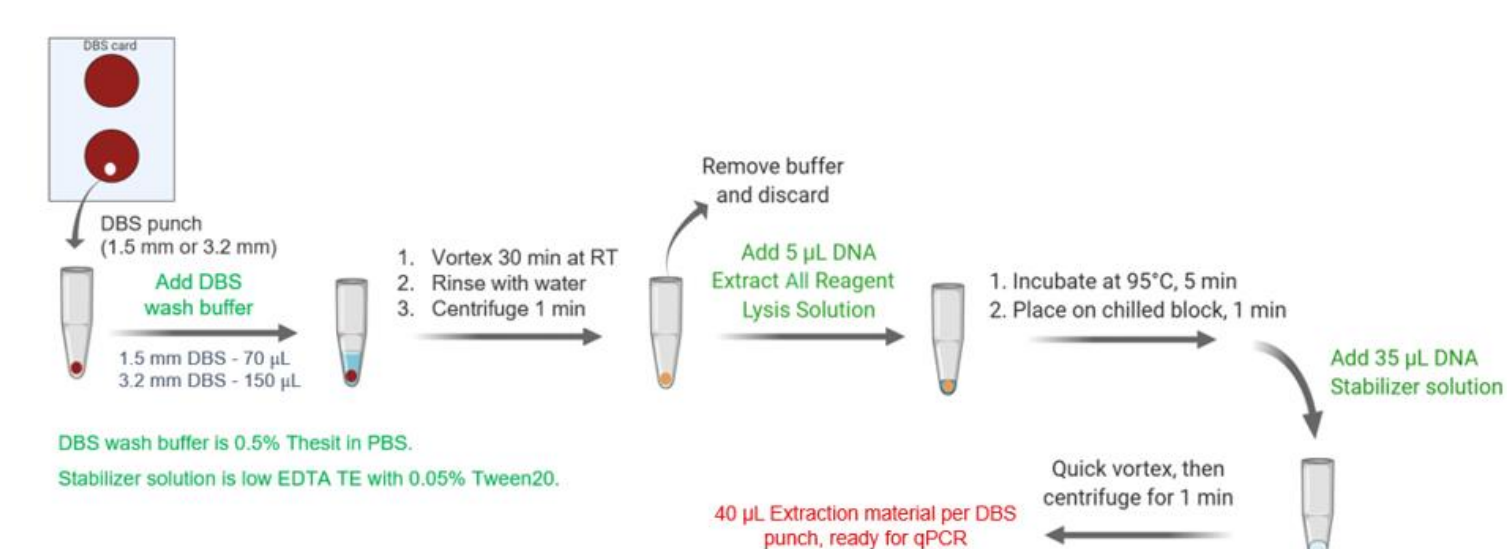
Intensive research in early detection of Severe Combined Immunodeficiencies (SCID) and Spinal Muscular Atrophy (SMA) is essential to discover prevention methods for infants' permanent disabilities or death. In addition to SCID, SMA has been recently added to the US-RUSP list. SMA is a motor neuron disorder caused by mutation in the SMN1 gene, whereas SCID constitutes a series of immune system functionality diseases exhibiting low levels of TREC and/or KREC. We have developed two versions of the TaqMan SCID/SMA multiplex assay (with and without KREC), that permit concomitant detection of both SMA and SCID targets, and each assay includes RNase P gene as an internal genomic control. Additionally, a simple and robust DBS sample preparation method was developed using DNA Extract All Lysis solution (Figure 1).

Multiplex TaqMan™ Assay Fluorescent Label Assignment:

TaqMan™ SCID/SMA assay (A47929): SMN1- FAM, TREC- JUN™, RNaseP (genomic control)- VIC™. Primers and TaqMan® dual labeled probes were synthesized by Thermo Fisher and formulated into a 20X TaqMan™ Assay.

TaqMan™ SCID/SMA Plus assay (A48569): SMN1- FAM, KREC- ABY™, TREC- JUN™, RNaseP (genomic control)- VIC™. Primers and TaqMan™ dual labeled probes were synthesized by Thermo Fisher and formulated into a 20X TaqMan™ Assay.

Figure 1



- Sample preparation workflow duration is ~40 min total with ~35 min hands-off time. When using freshly prepared, routine DBS samples as opposed to desiccated DBS specimens, sample preparation duration is ~20 minutes total.
- DBS punch-to-answer duration for 96-well plate format is < 2 hr.

RESULTS

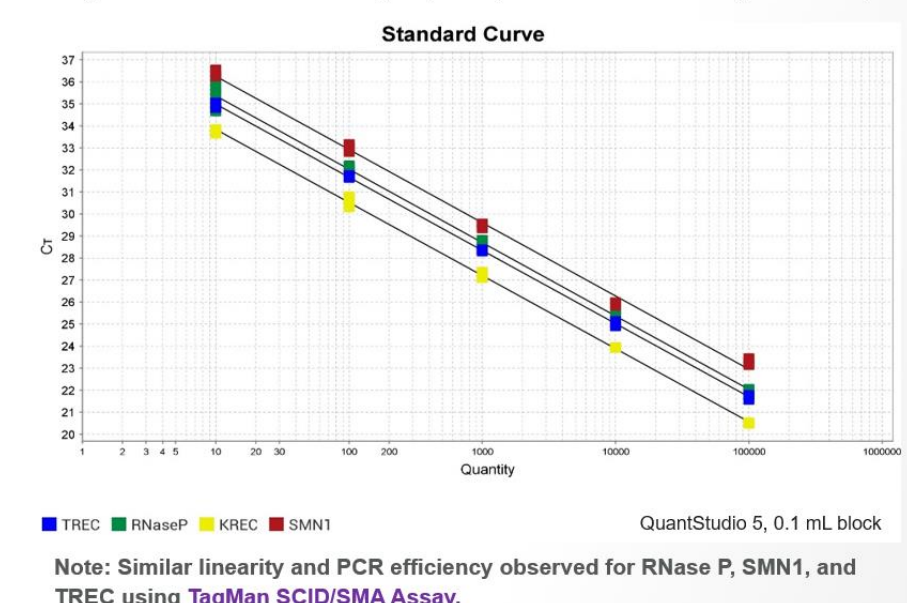
1. High linearity and PCR efficiency was observed for all targets (Figure 2). We demonstrate TREC and KREC copy number detection capability of as low as 10 copies per reaction and a LoQ of 25 copies per reaction.

Figure 2

Standard Curve

TaqMan SCID/SMA Plus Assay

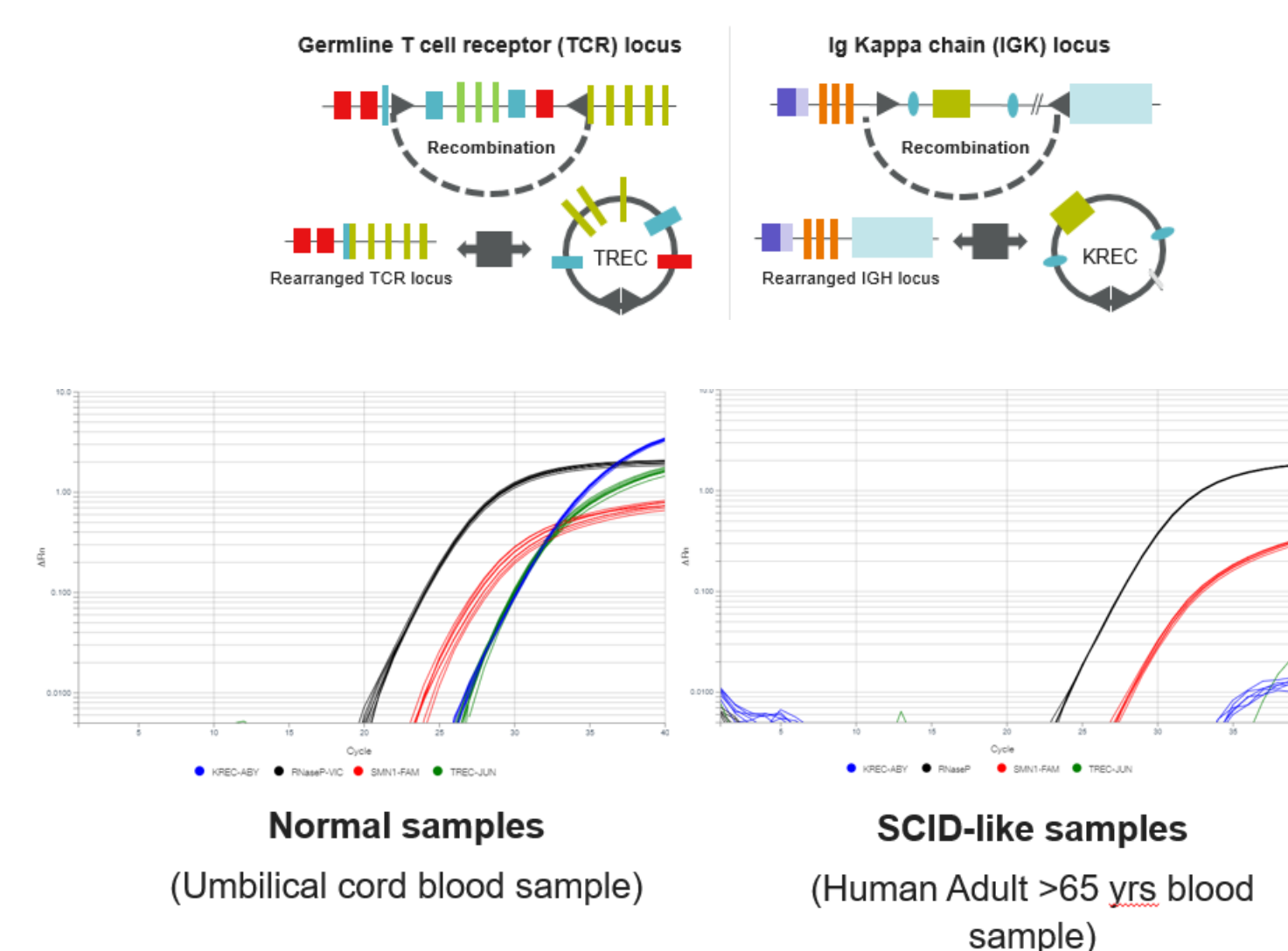
5 logs at 1:10 serial dilution of target specific plasmid construct starting at 1x10⁶ copies



LoD: 10 copies/reaction for RNase P, SMN1, TREC, KREC
LoQ: 25 copies/reaction for both TREC and KREC

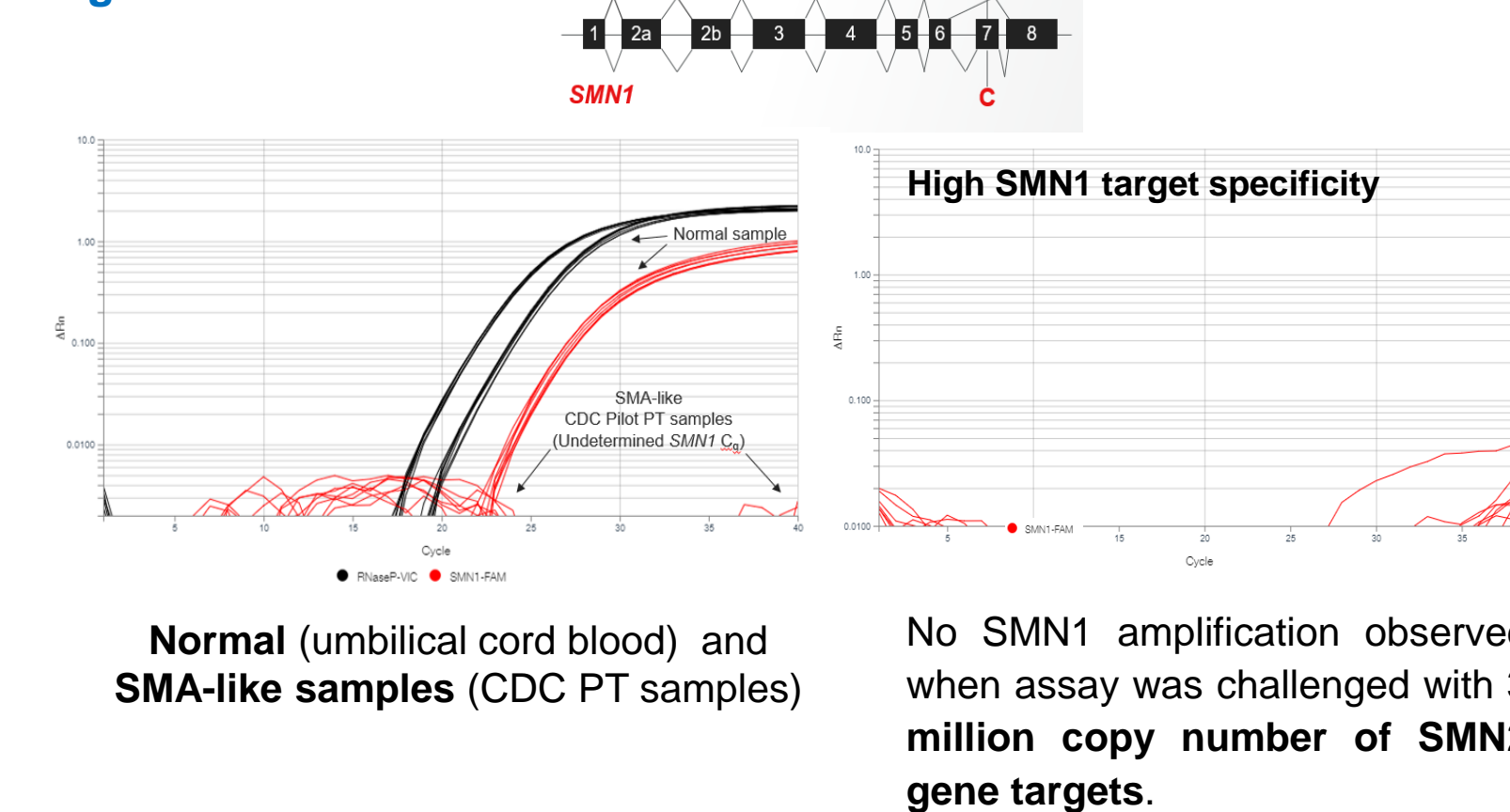
2. The TaqMan™ SCID/SMA Plus assay is designed to detect the presence of TREC and KREC (Figure 3), while the TaqMan™ SCID/SMA assay is designed to detect the presence of TREC only.

Figure 3



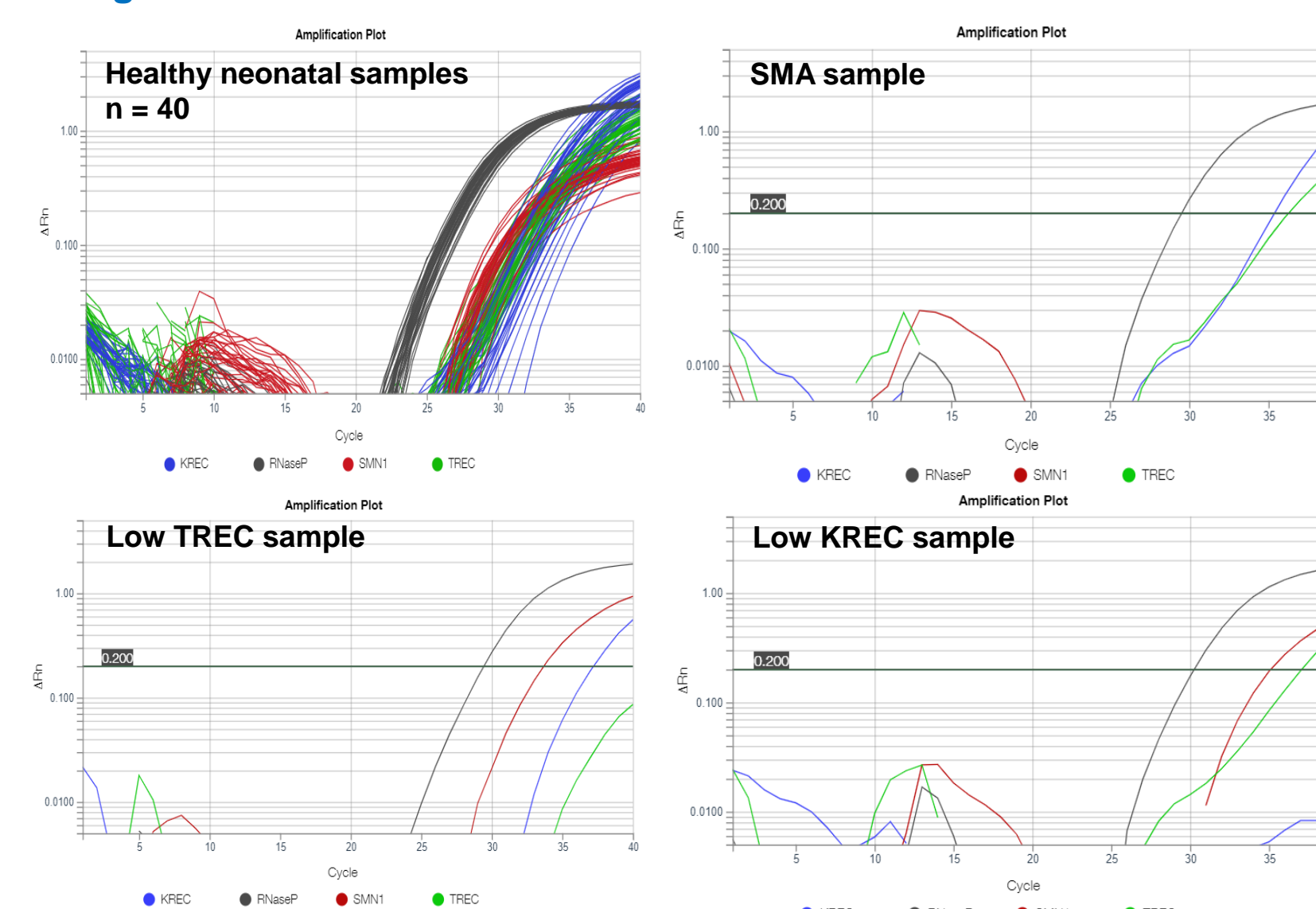
3. Both assays are designed to target exon 7 of SMN1 gene and effectively eliminated non-specific detection of the highly similar SMN2 gene (Figure 4 – generated with TaqMan™ SCID/SMA Plus assay).

Figure 4



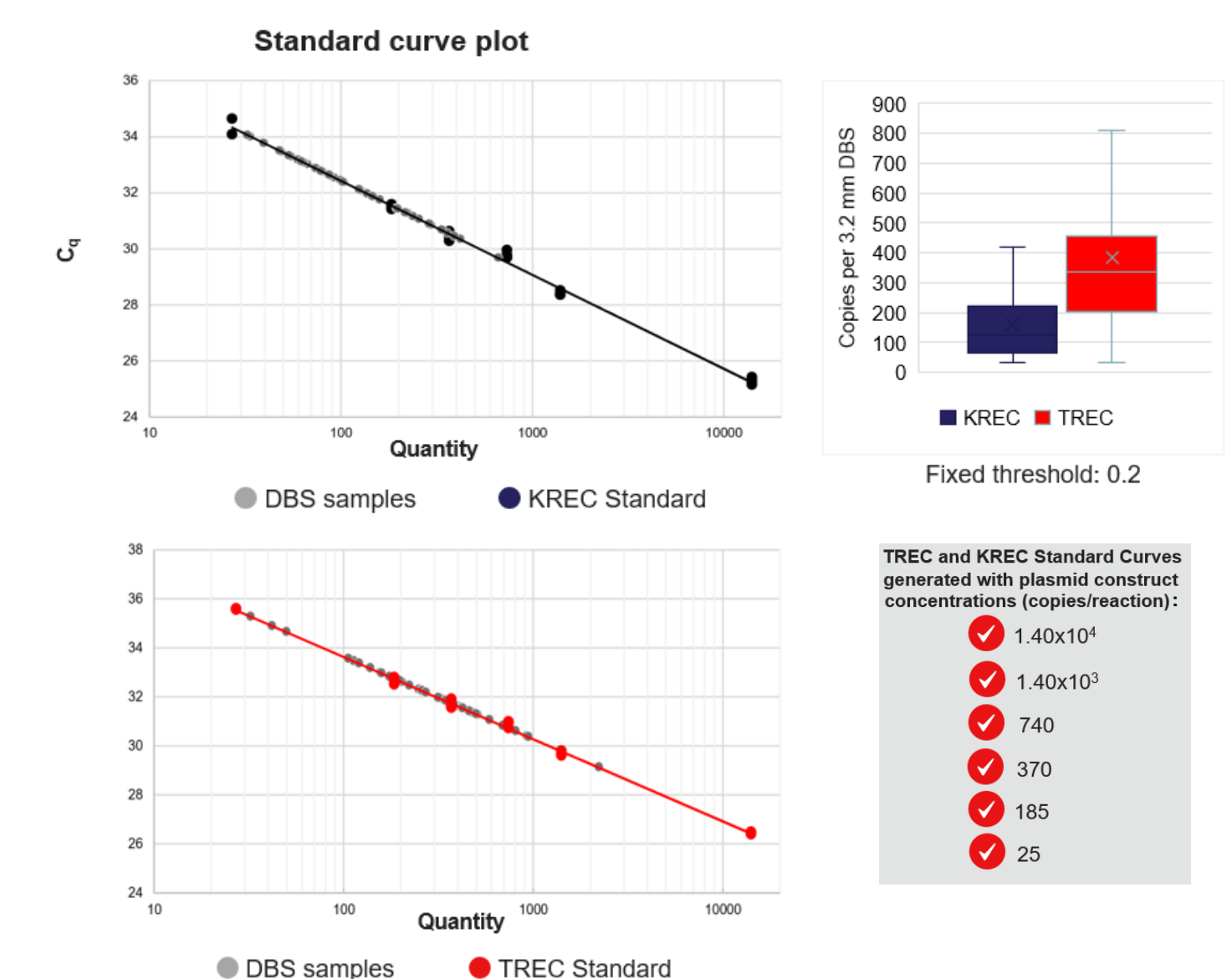
4. We demonstrate the performance of TaqMan™ SCID/SMA Plus assay with healthy neonatal DBS samples tested at Kumamoto University¹. Additionally, the assay's high analytical accuracy was evaluated using confirmed SMA positive samples and SCID positive samples containing low levels of TREC/KREC (Figure 5).

Figure 5



5. TREC and KREC quantitation: 46 clinical research DBS samples were quantified using the TaqMan™ SCID/SMA Plus Assay. TREC and KREC plasmid constructs were used to generate standard curves. See Figure 6.

Figure 6



CONCLUSION

In conclusion, we have developed a highly specific, sensitive, and robust multiplex assay for SMA and SCID testing with a rapid and streamlined turnaround workflow to aid further research efforts.

For newborn screening research



Recommended Applied Biosystems™ QuantStudio™ real-time PCR systems (96 and 384 blocks):

- QuantStudio™ 5 system
- QuantStudio™ 6 Flex system
- QuantStudio™ 7 Flex system
- QuantStudio™ Dx system (in RUO mode)
- QuantStudio™ 12K Flex system

Affiliations

1. Thermo Fisher Scientific, 3450 Central Express Way, Santa Clara, CA
2. Kumamoto University Graduate School of Medical Sciences, Department of Pediatrics

Author Contact Info

Sonu Baral, sonu.baral@thermofisher.com

Acknowledgements

We thank the CDC NBS team, Suzanne Cordovado et. al., for providing the Proficiency Testing (PT) blood card samples as well as for ongoing scientific discussion. We thank Prof. Nakamura, Dr. Sawada, and Kumamoto University lab team for providing data presented in this poster (see Figure 5). We also thank the Thermo Fisher Pleasanton oligo manufacturing for the primers and probes development and manufacturing.

Product links

TaqMan™ SCID/SMA assay:
<https://www.thermofisher.com/order/catalog/product/A47927#/A47927>

TaqMan™ SCID/SMA Plus assay:
<https://www.thermofisher.com/order/catalog/product/A48566#/A48566>