

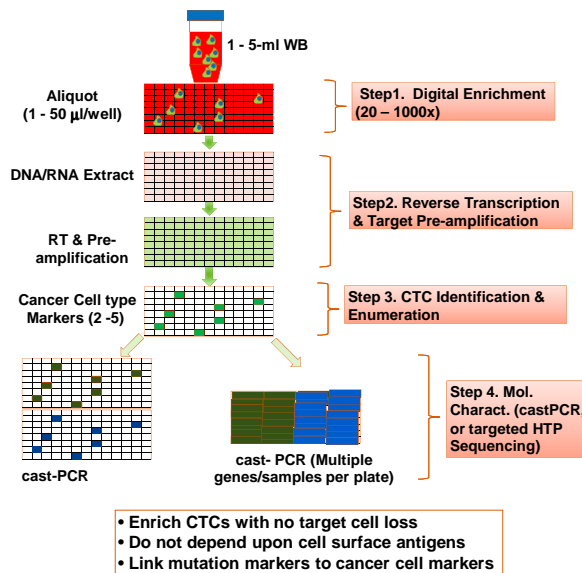
Digital Sample Enrichment and RT-castPCR Detection for Direct Molecular Characterization and Enumeration of Circulating Tumor Cells (CTC)

David Xingfei Deng, Scott Sproul, Yun Bao, Yu Wang, Fawn Wang, David Merrill, Pius Brzoska, and Caifu Chen
Life Technologies, 850 Lincoln Centre Dr. Foster City, CA94404, USA

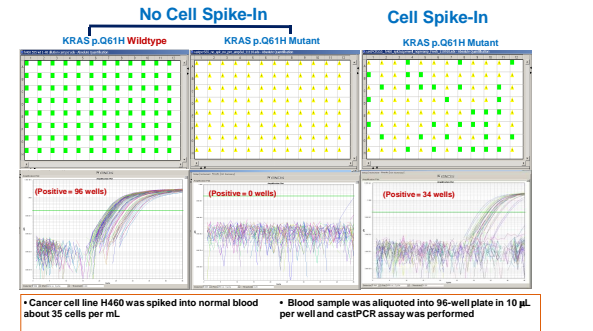


ABSTRACT Enumeration and molecular characterization of circulating tumor cells (CTCs) promise to be valuable for cancer diagnosis, survival prognosis, and treatment guidance. However, current methods require extensive enrichment process before analyzing rare CTCs in human blood. We reported here a new approach for direct CTC detection in whole blood samples by using digital sample enrichment and competitive allele-specific TaqMan PCR (castPCR) for rare mutations and RT-qPCR for cell type-specific genes. Whole blood samples from lung cancer patients or normal individuals with spiked-in known lung cancer cell lines were partitioned in aliquots of 2.5 - 50 μ L onto 96- or 384-well plates, such that each well contained either one cancer cell or none in the presence of 2×10^4 - 4×10^5 normal white blood cells and 10^7 - 2×10^8 red blood cells. The sample partition process resulted in a digital enrichment of 20 - 400 folds (the relative ratio of CTC to normal cells) in a CTC-positive well. castPCR clearly identified known mutations and CK19 in spiked-in samples of ~10 - 30 cells per mL whole blood, but there was no positive well in the absence of spiked-in cells. Furthermore, cell type specific markers (CK19) and known EGFR mutations were identified in the same sample wells, indicating that identified mutation was specifically derived from cancer cells. In two blood samples from lung cancer patients, EGFR mutation (p.L858R) was detected in all samples. Approximately, 50% of circulating lung tumor cells in a patient with positive EGFR p.L858R mutation had also positive EGFR p.T790M mutation, an inducible drug-resistant CTC marker. For those samples with negative detection of EGFR mutation, corresponding wild type sequences were detected in all sample wells. In conclusion, our data suggest that combination of digital sample enrichment with castPCR and RT-qPCR could be used to directly enumerate CTCs and detect cancer mutations in whole blood without prior biophysical sample enrichment. This new approach may pave the way for noninvasive CTC monitoring and individualized therapy.

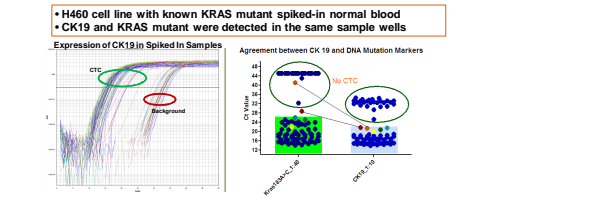
Digital Sample Enrichment and RT-castPCR Detection



castPCR Detection Accurately Determine Spiked-In Cancer Cells



Cell Type Markers (CK19) Co-relate To Mutation Markers



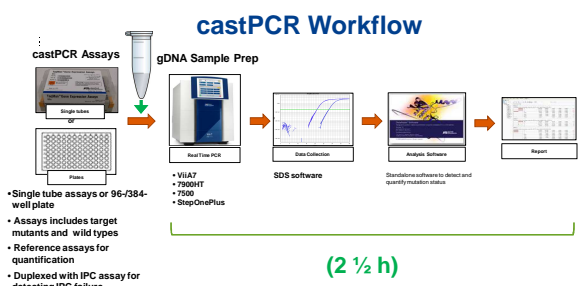
Challenges of CTC Analyses

- **CTCs Are Rare in Blood**
 - ~ 1 cell/mL Blood
 - Huge background cells (~1 in 10^6 WBC)
- **Variable Level of CTC Surface Antigen**
 - Unreliable isolation via cell surface markers
- **Ectopic Expression and Genetic Mutations in "Normal" Blood Cells**
 - Difficult in data interpretation from whole blood

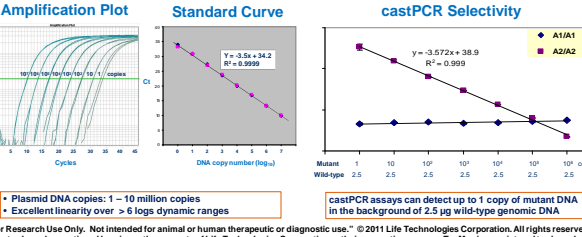
Our New Strategies

- **Sample Partition (Digital Enrichment)**
 - 100 - 1000 x enrichment
 - Independent of cell surface markers
- **High Selective Assays: castPCR**
 - Up to 1 in 10^6 selectivity
- **Mutation Detection from Both RNA and DNA**
 - Multiple copies in a cell
- **Co-relate Cancer cell markers and mutation markers in the same CTC**

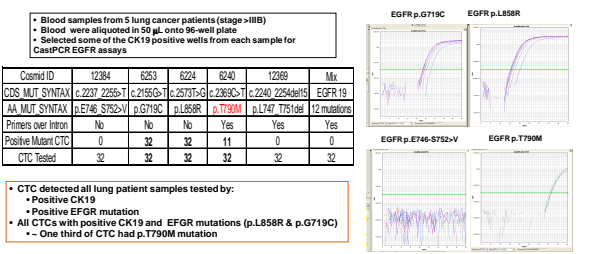
CastPCR Detection Has High Sensitivity and Specificity



castPCR Sensitivity and Selectivity



CastPCR Detection Accurately Determine CTCs From Lung Cancer



Summary

We present results for direct CTC molecular characterization and enumeration in blood:

- Up to 1000 fold sample enrichment by simple partition
- castRT-qPCR assays detecting cancer markers and mutations from the same CTCs
- New approach can be applied to CTCs of all cancer types for direct molecular characterization and enumeration

For Research Use Only. Not intended for animal or human therapeutic or diagnostic use. © 2011 Life Technologies Corporation. All rights reserved. The trademarks mentioned herein are the property of Life Technologies Corporation or their respective owners. TaqMan is a registered trademark of Roche Molecular Systems, Inc.