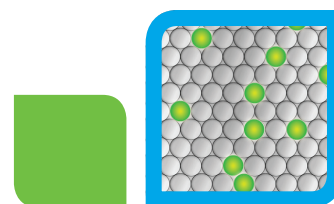


Droplet Digital™ PCR: TRAPing Telomerase Activity Using Droplet Digital PCR

Dawne N. Shelton,¹ Jack Regan,¹ George Karlin-Neumann,¹ Jue Lin,² Elizabeth Blackburn²
¹Digital Biology Center, Bio-Rad Laboratories, Inc., 5731 W Las Positas Blvd, Pleasanton, CA 94588
²Department of Biochemistry and Biophysics, UCSF School of Medicine, San Francisco, CA



Droplet Digital PCR

Bulletin 6541

Abstract

Telomerase is responsible for adding hexameric repeats to telomeres, which rewinds the mitotic clock and enables a cell to continuously divide. Abundant telomerase activity is found in fetal and adult stem cells, germ cells, and cancer. It is also present at much lower levels in nonpluripotent cells, such as immune cells, but these levels are difficult to measure using current methods.

The telomerase repeat amplification protocol (TRAP) measures the presence of active telomerase by measuring the activity of the enzyme on a starting template, which is then amplified by PCR. The current most sensitive method of detection still uses radioactivity and laborious PAGE sequencing gels followed by densitometry to quantify telomerase.

Here we explore using Droplet Digital PCR (ddPCR™) technology and EvaGreen chemistry to provide absolute quantification of telomerase activity. This was achieved through single-molecule counting of telomerase-extended templates that were partitioned into droplets, amplified by PCR, and detected by fluorescence droplet flow cytometry. Analysis of control samples suggests ddPCR is at least an order of magnitude more sensitive than TRAP radiography and is more amenable to higher-throughput analysis and lower input sample. This method extends throughput, sensitivity, and the range of biological samples that can be analyzed for telomerase activity.

Introduction

Telomeres are the protective structures at the ends of chromosomes and consist of 6 bp repeat sequences. In young cells, these regions can be as long as 15 kb and the telomeres act as caps, which protect the DNA ends. These ends naturally degrade with each passing cell division, usually losing 25–200 base pairs per division. Once they are shortened below a critical length (estimated to be 200–300 bp) the cells arrest and become senescent, or “old.” Telomeres can be thought of as a cellular or mitotic clock. Once the clock has wound down, the cells either die or pass through crisis and become immortal.

One of the mechanisms of immortality is activation of the enzyme telomerase. Telomerase is a reverse transcriptase enzyme that adds hexameric repeats to the ends of telomeres using an internal RNA template. Extension of telomeres allows

cells to proliferate. The vast majority of adult cells do not express telomerase. Adult stem cells, germ cells, and cancer cells are the exceptions.

The aim of this work was to develop a more sensitive high-throughput assay for measuring telomerase activity.

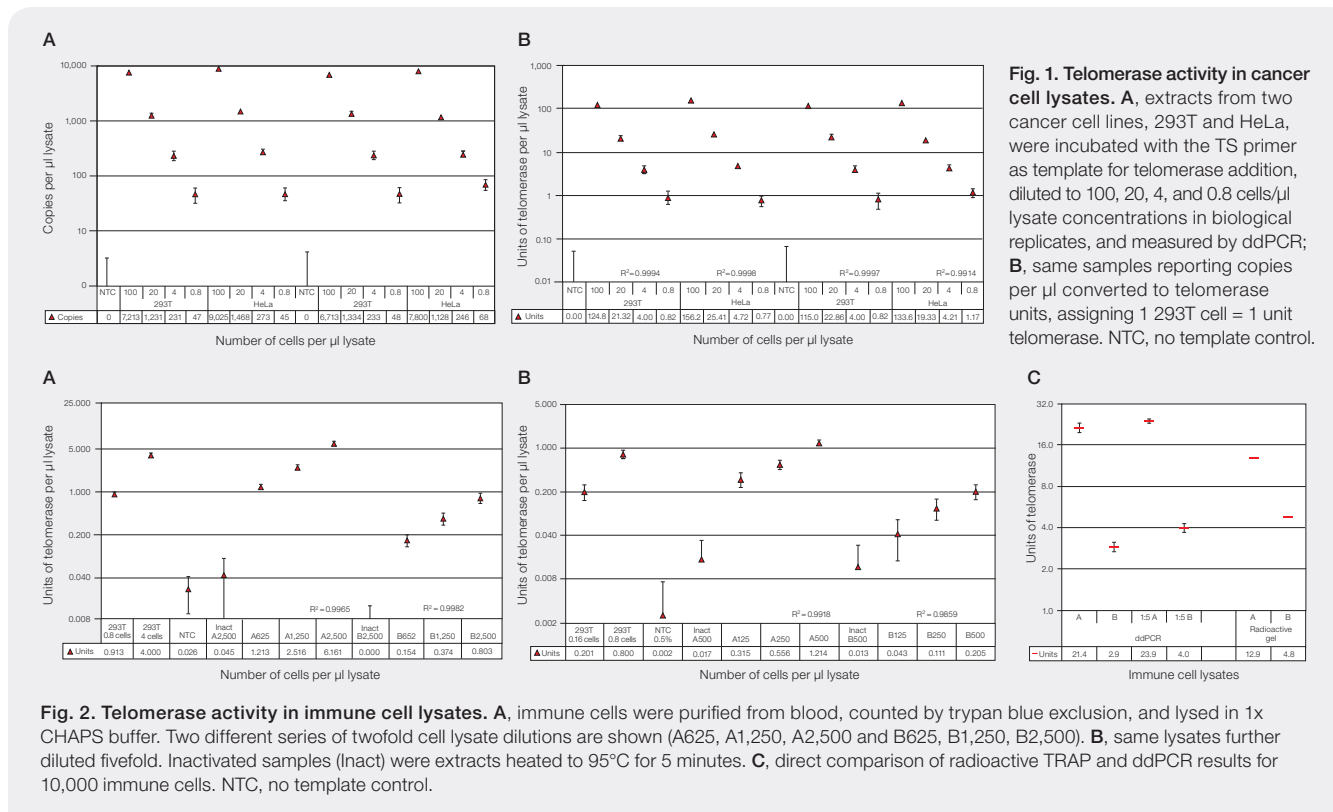
Methods and Results

Complementary DNA was made from cell lysates containing active telomerase enzyme and a starting template, the TS oligonucleotide. Telomerase-extended templates were partitioned into droplets, amplified by PCR, and detected by fluorescence droplet flow cytometry using the QX200™ Droplet Digital PCR system and EvaGreen. We investigated two commonly used cancer cell lines, HeLa and 293T, for telomerase activity in parallel fivefold dilution series in duplicate. Templates were counted by ddPCR, and units of telomerase were calculated based on 1 293T cell = 1 unit of

telomerase (Figure 1). Quantities of telomerase ranged from 0.77 units in 0.8 cells per reaction to 156 units in 100 cells per reaction. R² values for each series ranged from 0.9914 to 0.9998, demonstrating excellent linearity.

To further test the range of the system, lysates from purified normal immune cells were examined in two sets of twofold dilution series from two patient samples. Again linearity and

separation from controls were excellent. R² values for each series ranged from 0.9859 at the lowest levels to 0.9982 at the higher telomerase levels. Patient sample B expressed 3–4 units of telomerase per 10,000 cells while patient A exhibited 21–24 units per 10,000 cells. We were able to clearly detect 0.043 units of telomerase in 125 cells of patient B, demonstrating at least 5 orders of sensitivity (Figure 2).



Conclusions

- Droplet Digital PCR provides precise and accurate concentration measurements of telomerase activity on an absolute scale
- Analysis of control samples suggests ddPCR is more sensitive and linear than TRAP radiography
- This method extends throughput, sensitivity, and the range of biological samples that can be analyzed for telomerase activity

For more information, visit

www.bio-rad.com/web/ddPCRTRAP.

EvaGreen is a trademark of Biotium, Inc.

The QX200 Droplet Digital PCR system and/or its use is covered by claims of U.S. patents, and/or pending U.S. and non-U.S. patent applications owned by or under license to Bio-Rad Laboratories, Inc. Purchase of the product includes a limited, non-transferable right under such intellectual property for use of the product for internal research purposes only. No rights are granted for diagnostic uses. No rights are granted for use of the product for commercial applications of any kind, including but not limited to manufacturing, quality control, or commercial services, such as contract services or fee for services. Information concerning a license for such uses can be obtained from Bio-Rad Laboratories. It is the responsibility of the purchaser/end user to acquire any additional intellectual property rights that may be required.



Bio-Rad
Laboratories, Inc.

Life Science
Group

Web site www.bio-rad.com USA 800 424 6723 Australia 61 2 9914 2800 Austria 01 877 89 01 Belgium 09 385 55 11 Brazil 55 11 3065 7550
Canada 905 364 3435 China 86 21 6169 8500 Czech Republic 420 241 430 532 Denmark 44 52 10 00 Finland 09 804 22 00
France 01 47 95 69 65 Germany 089 31 884 0 Greece 30 210 9532 220 Hong Kong 852 2789 3300 Hungary 36 1 459 6100 India 91 124 4029300
Israel 03 963 6050 Italy 39 02 216091 Japan 81 3 6361 7000 Korea 82 2 3473 4460 Mexico 52 555 488 7670 The Netherlands 0318 540666
New Zealand 64 9 415 2280 Norway 23 38 41 30 Poland 48 22 331 99 99 Portugal 351 21 472 7700 Russia 7 495 721 14 04
Singapore 65 6415 3188 South Africa 27 861 246 723 Spain 34 91 590 5200 Sweden 08 555 12700 Switzerland 026674 55 05
Taiwan 886 2 2578 7189 Thailand 1800 88 22 88 United Kingdom 020 8328 2000