

SuperScript® IV Reverse Transcriptase: A New Reverse Transcriptase for RNA Analysis

invitrogen
by Thermo Fisher Scientific

Authors: Bianca J Lam¹, Kevin Zhang¹, Joanna Guo¹, Lushen Li¹, Kimberly Wong Garcia¹, Asta Jokubauskaite², Milda Kaniusaite², Tomas Radzvilavicius², and Arunas Lagunavicius²

¹Thermo Fisher Scientific/Life Sciences Solutions, 5781 Van Allen Way, Carlsbad, California, USA, 92008

²Thermo Fisher Scientific/Life Sciences Solutions, V.A. Graiciuno 8, LT-02241 Vilnius, Lithuania

ABSTRACT

Survey and interview studies conducted over a three year period revealed that researchers are not satisfied with their current reverse transcriptase and are performing reactions with increasingly difficult samples, such as poorly purified RNA and unpurified RNA (direct RT) that both contain inhibitors. To meet this performance gap, the Thermo Fisher Life Sciences Solutions group produced a new reverse transcriptase, SuperScript® IV, and experiments we performed show that it is the most robust reverse transcriptase compared to other enzymes. SuperScript® IV characterization was performed in the context of "real world" situations where users do not have perfect RNA samples. In the presence of a variety of inhibitors, we demonstrate that SuperScript® IV possesses superior performance in a variety of inhibitors, such as alcohols, salts, detergents, phenol, heparin, hematin, bile salts, and formalin typically found in sample preparation reagents, cell lines, blood, feces, and FFPE samples. This enzyme can even detect RNA targets in unpurified RNA samples (directly lysed cells) and whole blood without sacrificing sensitivity and yield. The introduction of SuperScript® IV enables researchers to obtain more consistent results independent of sample quality and simplify and speed up workflows by eliminating RNA purification.

INTRODUCTION

Gene expression starts when RNA is transcribed from DNA. Expression levels of different RNA targets can be used to characterize species, tissue types, cell types, and healthy and diseased cells. Because RNA is unstable, it is necessary to convert these molecules into more stable ones without loss of quantitative and coding information. Thus, reverse transcriptases are indispensable enzymes because they convert RNA into DNA (cDNA), a stable form of the genetic code. SuperScript® IV is the newest member of the SuperScript® family of reverse transcriptases, which is known for its quality and reliability in cDNA synthesis. This new enzyme has been proven to function in experimentally challenging conditions that many scientists face, such as in the presence of enzyme inhibitors found in reagents, cells, blood, feces, and FFPE samples. To demonstrate this enzyme's capabilities even further, RNA transcripts were detected in unpurified RNA samples.

REVERSE TRANSCRIPTION PROTOCOL

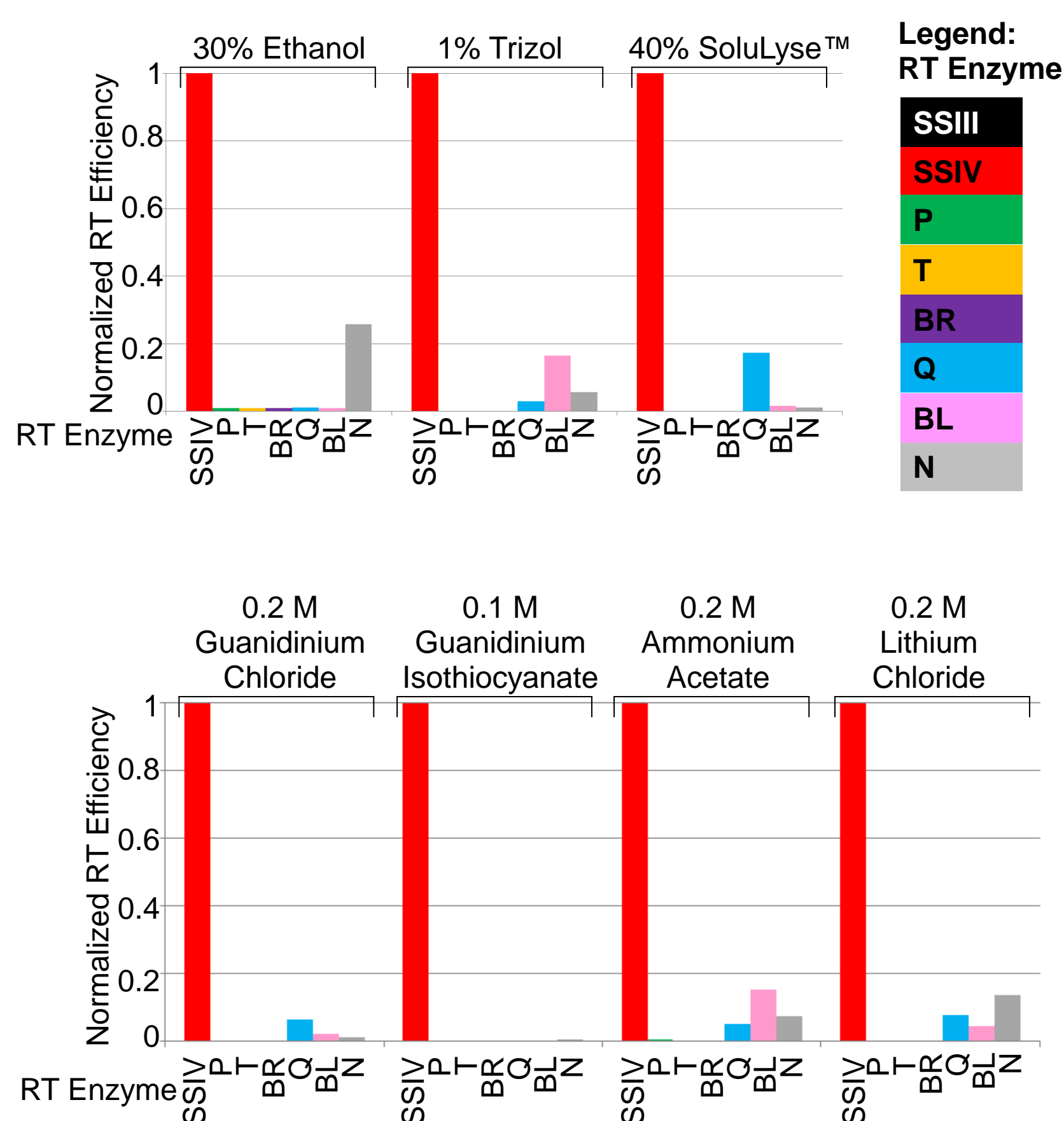
Step	Procedure	Procedure details
1	Annual primer to template RNA	<p>a. Combine the following components in a reaction tube:</p> <p>Note: Consider the volume for all components listed in steps 1 and 2 to determine the correct amount of water required to reach your final reaction volume.</p> <p>Component</p> <p>50 µM oligo-dT₁₈ primer, 50 µM random hexamers, or 2 µM gene-specific reverse primer 1 µL</p> <p>10 mM dNTP mix (10 mM each) 1 µL</p> <p>Template RNA (10 µg total RNA or 10 µg-500 ng mRNA) up to 11 µL</p> <p>100 mM RNase inhibitor 1 µL</p> <p>100 mM DTT 1 µL</p> <p>100 mM EDTA 1 µL</p> <p>100 mM Tris-HCl, pH 8.0 1 µL</p> <p>b. Mix and briefly centrifuge the components.</p> <p>c. Heat the RNA-primase mix at 65°C for 5 minutes, and then incubate on ice for at least 1 minute.</p>
2	Prepare RT reaction mix	<p>a. Mix and briefly centrifuge the 5x RT Buffer.</p> <p>b. Combine the following components in a reaction tube:</p> <p>Component</p> <p>5x RT Buffer 4 µL</p> <p>100 mM DTT 1 µL</p> <p>100 mM EDTA 1 µL</p> <p>100 mM Tris-HCl, pH 8.0 1 µL</p> <p>SuperScript® IV Reverse Transcriptase (200 U/µL) 1 µL</p> <p>c. Cap the tube, mix, and then briefly centrifuge the contents.</p>
3	Combine annealed RNA and RT reaction mix	<p>Add RT reaction mix to the annealed RNA.</p>
4	Incubate reactions	<p>a. If using random hexamers, incubate the combined reaction mixture at 25°C for 10 minutes, and then proceed to step b.</p> <p>b. If using oligo-dT₁₈ or gene-specific primers, directly proceed to step b.</p> <p>c. Incubate the combined reaction mixture at 50-55°C for 10 minutes.</p> <p>d. Inactivate the reaction by incubating at 85°C for 10 minutes.</p>
5	Optional: Remove RNA	<p>Note: Amplification of some PCR targets (2-3 kb) may require removal of RNA. To remove RNA, add 1 µL of 10% SDS and incubate at 37°C for 20 minutes.</p>
6	PCR amplification	<p>Use your RT reaction immediately for PCR amplification or store it at -20°C.</p> <p>Note: As a recommended starting point for PCR, reverse transcription reaction (cDNA) should contain 10% of the total reaction volume.</p>

RESULTS

Figure 1. Reverse transcriptase inhibitors

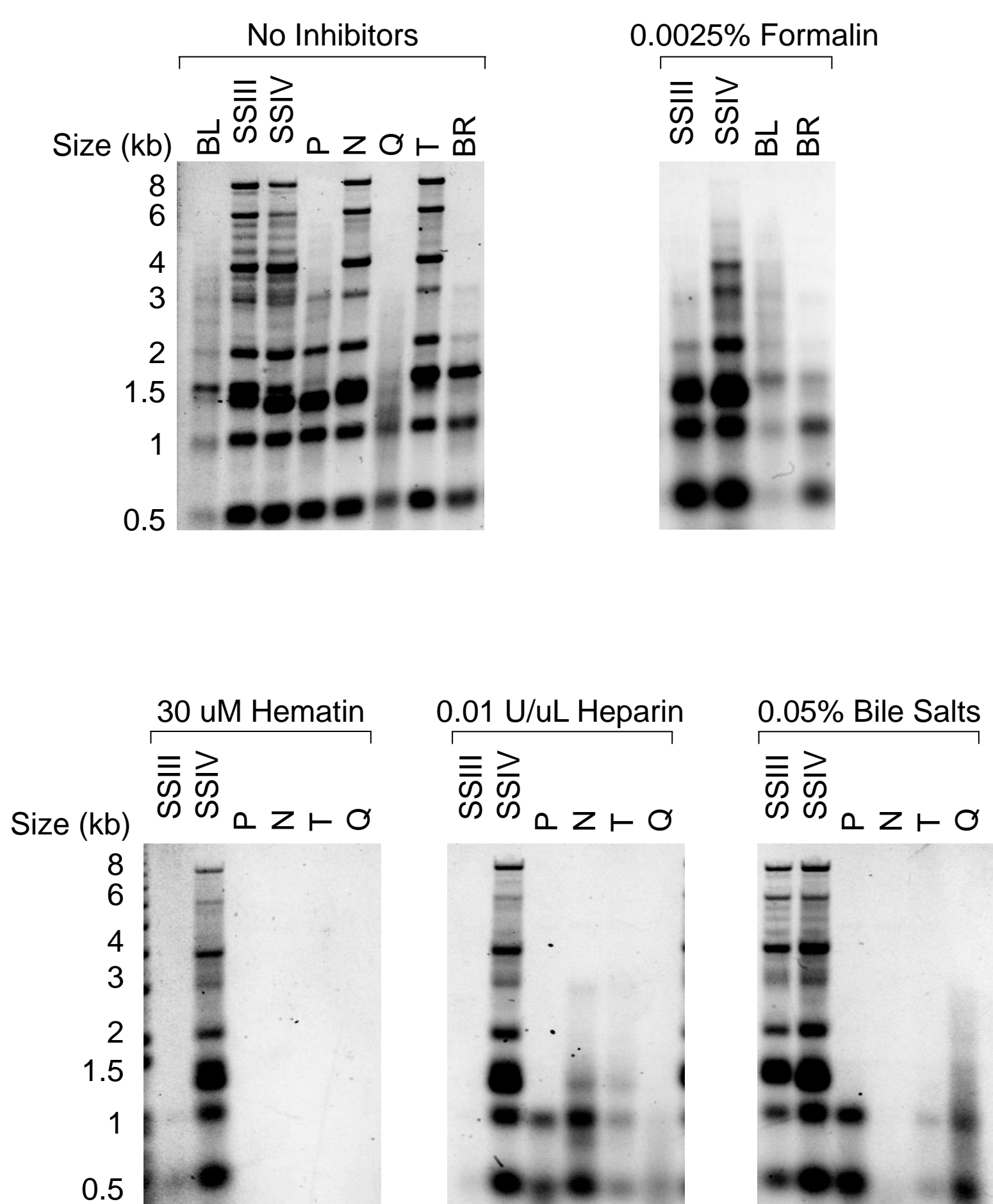
Inhibitors	Source
Alcohols (Ethanol)	Sample prep
Salts	Sample prep
• Guanidinium chloride	
• Guanidinium isothiocyanate	
• Ammonium acetate	
• Lithium chloride	
Cell lysis	Sample prep
• Trizol	
• SoluLyse	
Heparin, hematin, bile salts	Cells, blood, feces
Formalin	FFPE

Figure 2. SuperScript® IV RT-qPCR Performance in the Presence of Inhibitors



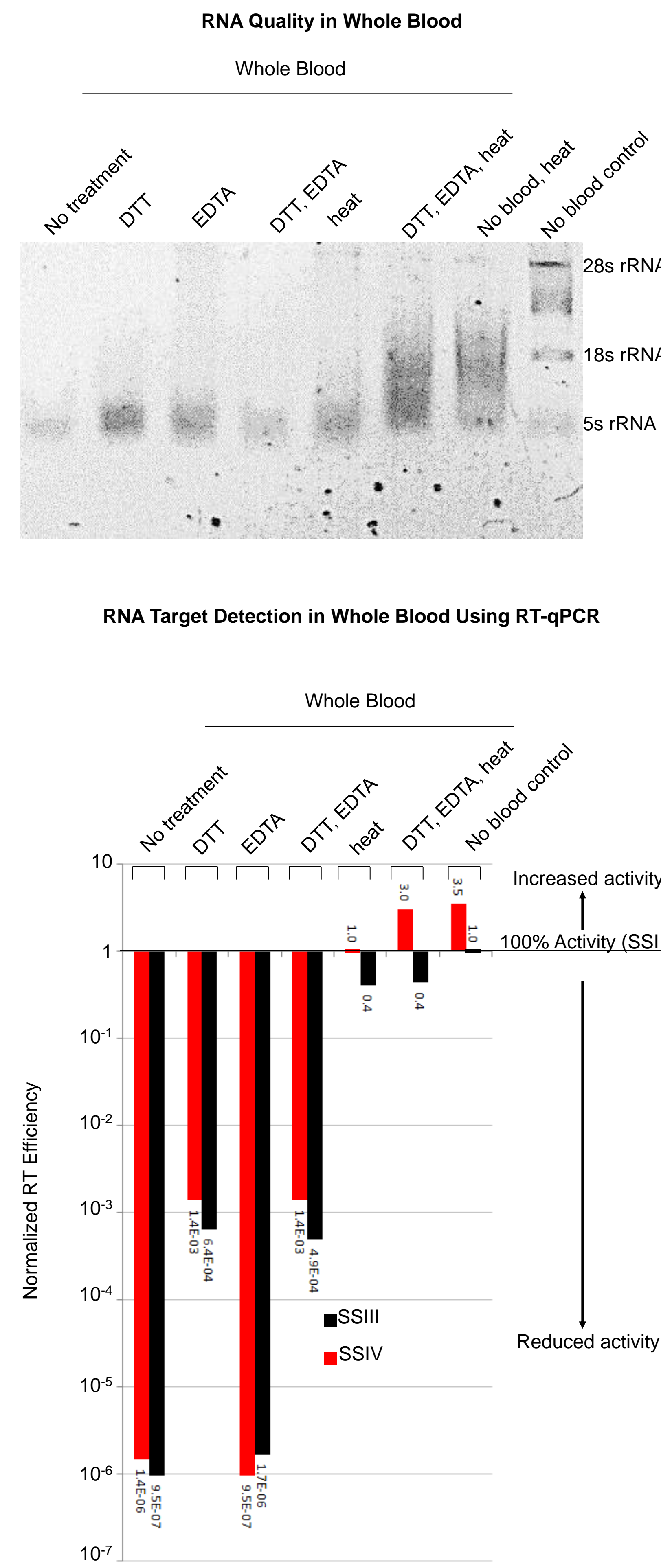
100 ng of total RNA from HeLa cells (Cat# AM7852) was used in a 20-µL SSIV reaction with oligo(dT)₂₀ according to provided protocol. Competitive products followed the manufacturer's recommended protocols. Inhibitors were added to total RNA prior to annealing to primer or addition of RT reaction mix. 10% of cDNA in the RT reaction composed 10% of the TaqMan® qPCR (Cat# 11785-01K) reaction volume which targeted the human TFRC (CD71) (transferrin receptor) gene (Cat# 4333770F). Ct values were normalized to SuperScript® IV using the equation: Normalized Y values = $2^{-(Ct_{SSIV} - Ct_{SSIII})}$. SSIV consistently detects targets in HeLa RNA in the presence of inhibitors commonly found in sample preparation reagents and blood.

Figure 3. SuperScript® IV First-Strand cDNA Performance in the Presence of Inhibitors



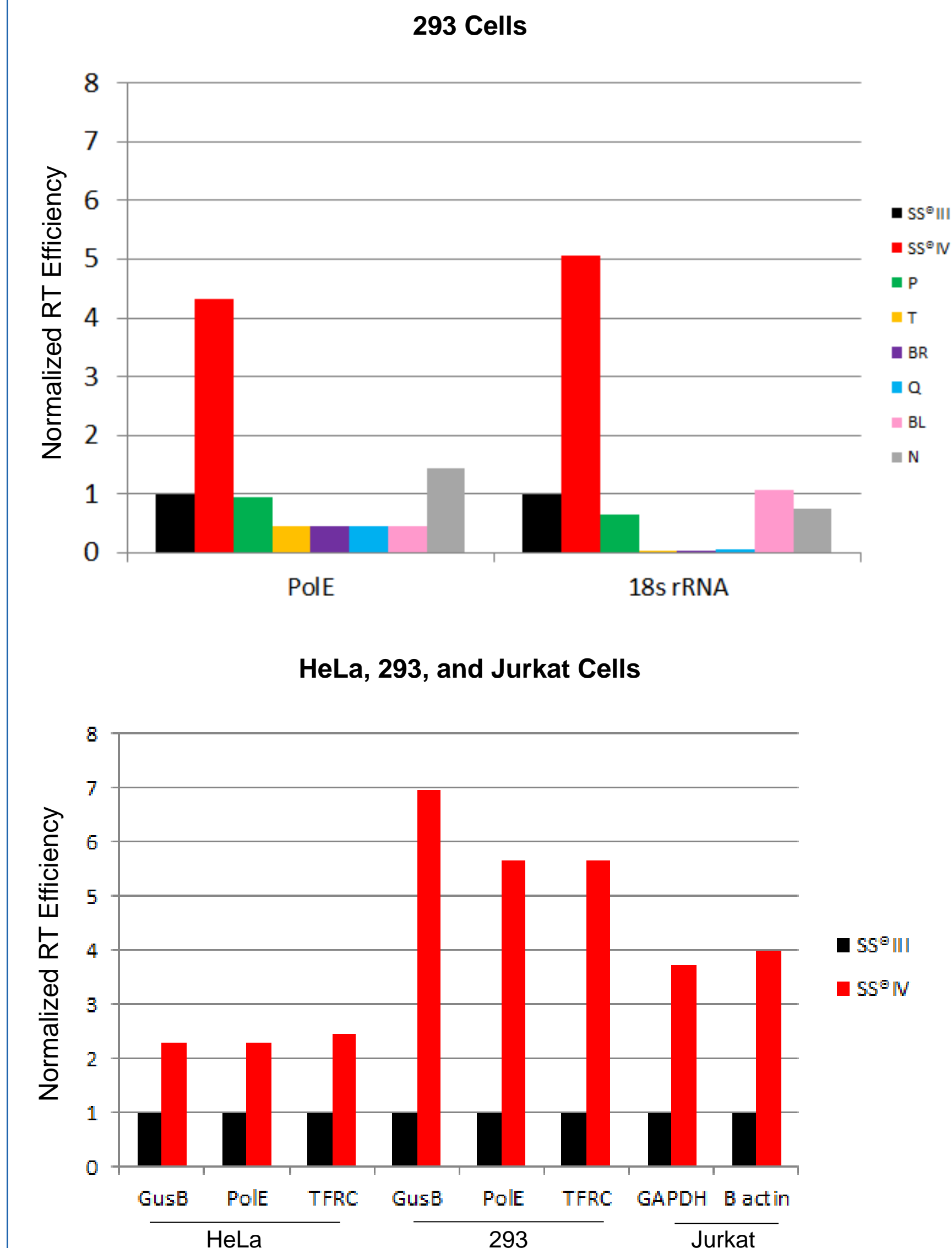
500 ng 0.5-10 kb RNA ladder (Cat# 15623-200) was used in a 10-µL SSIV reaction with oligo(dT)₂₀ according to provided protocol. Competitive products followed the manufacturer's recommended protocols. Inhibitors were added to total RNA prior to annealing to primer or addition of RT reaction mix. First-strand cDNA were resolved by alkaline gel electrophoresis and cDNA was stained using SYBR-gold (Cat# S-11494). NaOH hydrolyzes all RNA resulting in only visualization of cDNA. SSIV consistently generates longer cDNA with higher yields in the presence of inhibitors found in FFPE (formalin), blood (hematin and heparin), and feces (bile salts).

Figure 4. RNA Detection in Whole Blood using SuperScript® IV



Human disease research may sometimes require the detection of RNA in whole blood. Whole blood contains a combination of inhibitors, including heparin, heme, and high concentrations of RNases. To investigate, the performance of SSIV in whole blood, 100 ng HeLa RNA was added to whole blood in reverse transcription reactions composed of 2.5% whole blood. Either untreated blood, addition of 5 mM DTT and/or 1 mM EDTA, and/or heat treatment was also tested to optimize RNA detection in blood. Heat treatment consisted of heating blood/RNA samples to 95°C for 10 minutes. As a positive control, the same amount of HeLa RNA was used in a reverse transcription reaction in the absence of blood. Normalized Y values = $2^{-(Ct_{SSIII} - Ct_{SSIV})}$. SSIV consistently detected more 18s rRNA target by TaqMan® qPCR in heat-treated whole blood than SSIII. The best results consisted of the addition of DTT, EDTA, heat treatment, and the use of SSIV, suggesting that the most inhibitory component in blood are RNases. Ct value with no blood is 14.

Figure 5. RNA Detection in Unpurified RNA



Pelleted cells were ground to a fine powder in liquid nitrogen. The powder was transferred to a microfuge tube and TE was added, vortexed, and centrifuged to pellet debris. Resulting clarified supernatant was transferred into a fresh microfuge tube. Prior to reverse transcription, EDTA and DTT were added to the supernatant to a final concentration of 1 mM and 5 mM respectively, and heated to 95 °C for 10 minutes. TaqMan® qPCR reactions were used to detect targets indicated in figures above. Ct values were normalized to SuperScript® IV using the equation: Normalized Y values = $2^{-(Ct_{SSIV} - Ct_{SSIII})}$. SSIV consistently detects targets in unpurified RNA samples better than other commercial reverse transcriptases. Ct values for SSIV ranged between 25-35.

CONCLUSIONS

Thermo Fisher Scientific enables researchers to rapidly advance their studies by continuously evolving its reagents and tools for current and future research. The SuperScript® IV reverse transcriptase is an example of such a reagent. This enzyme is efficient in detecting RNA from samples containing inhibitors, such as alcohols, cell lysis reagents, salts, and those inherent in biological samples, such as those found in blood, feces, and FFPE RNA. Not only is this enzyme resistant to inhibitors, but it also possesses increased sensitivity such that RNA targets are detected from directly lysed human cell lines. Therefore, SuperScript® IV provides the most sensitive and reliable RNA analyses from the most challenging types of sample sources. Future studies will focus on more systematic and quantitative studies using unpurified RNA. Future considerations relating to SuperScript® IV reverse transcriptase will investigate more convenient formats for users running high throughput RT-qPCR and RT-PCR applications.

ACKNOWLEDGEMENTS

We would like to thank Toby Jordan and Samantha Li for their work and support in the development of SuperScript® IV.

TRADEMARKS/LICENSING

© 2015 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified. TaqMan is a registered trademark of Roche Molecular Systems, Inc., used under permission and license. TRIzol is a trademark of Molecular Research Center, Inc. SoluLyse is a trademark of Genlantis, Inc.

Let no sample go **unconquered** Introducing all-new SuperScript® IV RT

