

# Quantification of 100 RNA Targets from Limited Samples and Single Cells

Steven Okino, Michelle Kong, and Yan Wang  
Bio-Rad Laboratories, Inc., 2000 Alfred Nobel Drive, Hercules, CA, 94547, USA

BIO-RAD

## Introduction

Gene expression profiling is often limited by small sample size and low target gene expression. To address these issues we developed SsoAdvanced™ PreAmplification Supermix (172-5160, Bio-Rad Laboratories), a PCR-based reagent that increases the cDNA levels of up to 100 target genes at least 1000-fold. To validate our reagent, we analyzed a panel of genes that are differentially regulated during stem cell differentiation. We find that relative gene expression results generated using this product are statistically equivalent to results obtained using standard qPCR; however, 10,000-fold less starting material can be used. Importantly, SsoAdvanced PreAmplification Supermix maintains patterns of gene expression changes across samples; thus the same biological insights would be derived from a pre-amplification (PreAmplification) experiment and a standard gene expression profiling experiment. In addition, we demonstrate the application of this product for analyzing gene expression in single cells.

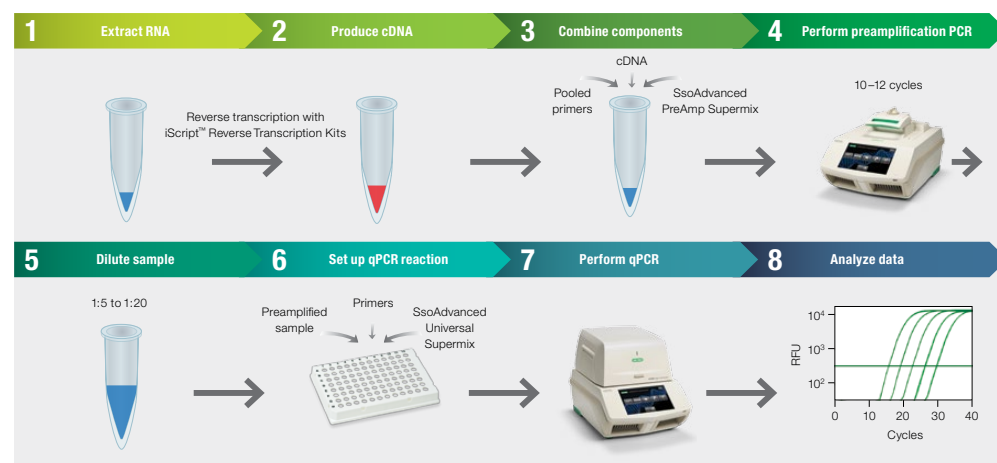
## Background

### Genes Analyzed

We analyzed a panel of 100 genes that are involved in stem cell differentiation. PrimePCR™ PreAmplification and qPCR assays were used in the PreAmplification reaction and for qPCR analysis, respectively.

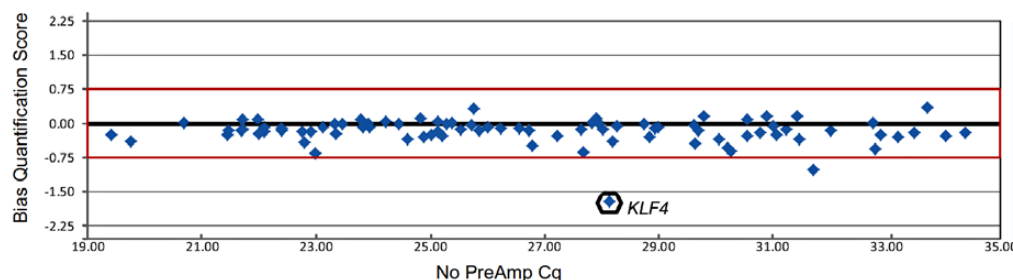
gDNA	POU5F1	ZFP42	CD44	DPPA5	GATA2	KLF4	NKX2-5	SMAD1	TAL1
HPRT1	POU3F2	DNMT3B	CDC42	EN2	GATA6	LIN28A	NROB1	SMAD2	TAT
B2M	NEUROD1	NRS2A2	CDK1	ENG	GDF3	MEIS1	NTSE	SMAD3	TCF3
GA PDH	LEFTY2	NR6A1	CHD1	ESRRB	GFAP	MESP1	OLIG2	SOX15	TCL1A
TFRC	UTF1	SOX2	CHD7	ETV2	GSC	MIXL1	OTX2	SOX17	TEK
TBP	TGDF1	ACTA2	CNOT3	FGF2	HAND1	MYBL2	PAF1	SOX3	TERT
NANOG	NEUROG2	AICDA	DES	FGF5	HNF4A	MYC	RIF1	SOX7	THAP11
PAX6	ASCL1	ALB	DPPA2	FLI1	HSPA9	MYOD1	ALA S1	STAT3	THY1
TBX3	ASCL2	ALP L	DPPA3	FOXP1	ISL1	NCAM1	SALL4	TBP	TRIM28
PRDM14	KLF2	CCNA2	DPPA4	GATA1	KAT5	NE S	SCN1A	TAGLN	ZFX

### Pre-amplification workflow



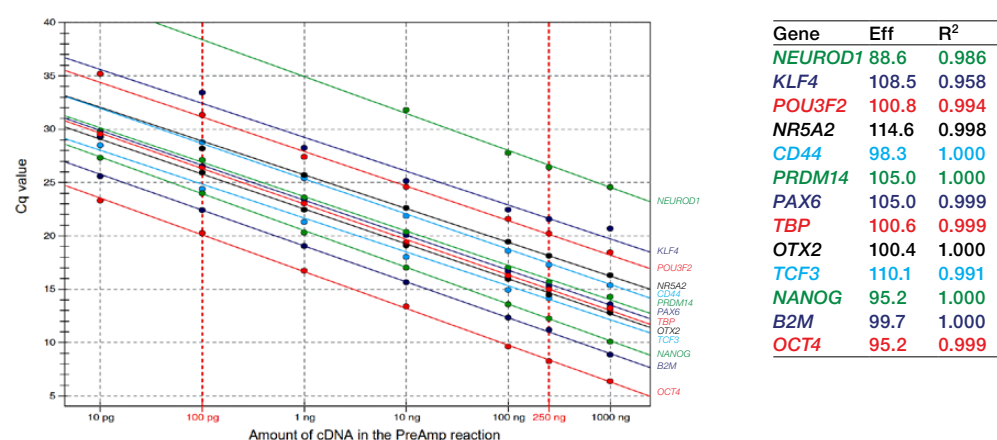
### SsoAdvanced PreAmplification Supermix generates minimal bias

To assess bias introduced by PreAmplification, we compared target gene levels in a cDNA sample before and after PreAmplification. For each target, we calculated the bias introduced by PreAmplification and plotted it against the no PreAmplification Cq values for the target genes. For 98% of genes with a no PreAmplification Cq < 35, PreAmplification did not introduce bias by more than 0.75 Cq.



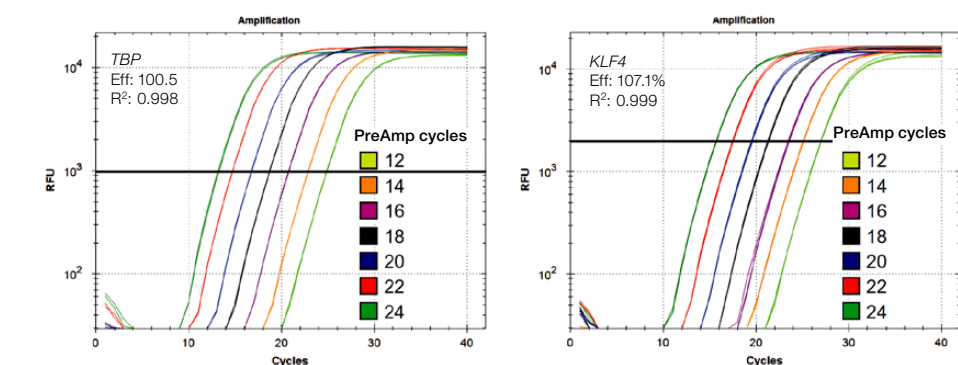
### SsoAdvanced PreAmplification Supermix has a wide dynamic range

PreAmplification was performed on 10 pg to 1 µg cDNA. Selected target genes with varied expression levels were analyzed by qPCR. The results (efficiency, R<sup>2</sup>) show that a wide linear dynamic range of cDNA input can be achieved regardless of target expression level.



### SsoAdvanced PreAmplification Supermix is efficient for up to 24 PreAmplification cycles

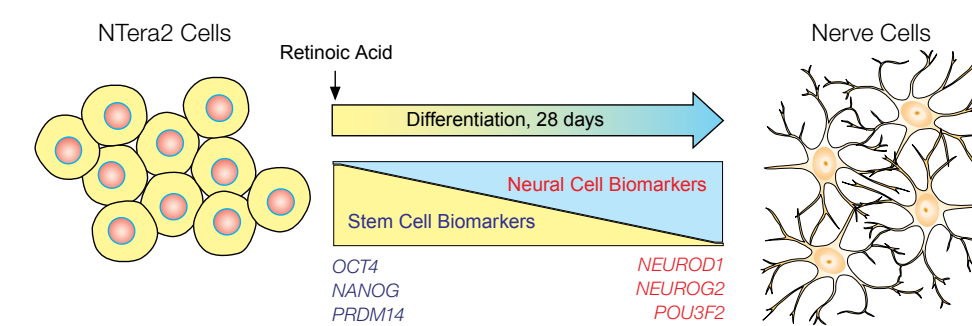
PreAmplification was performed on 10 ng cDNA for 12-24 cycles. qPCR traces for TBP and KLF4 are shown. Efficiency and R<sup>2</sup> values show that this product maintains exponential amplification for up to 24 PreAmplification cycles.



## Gene expression profiling during differentiation

### Analysis of cell populations undergoing differentiation

NTera2 cells (NT2) are a well-established model system used to study human stem cell behavior (Andrews 1984). When treated with retinoic acid (RA), NT2 differentiate into neurons (Pleasure and Lee 1993). As NT2 differentiate, the level of stem cell specific biomarkers decrease and the level of neuron-specific biomarkers increase (Deb-Rinker et al. 2005). We used the NT2 model system to determine if PreAmplification can accurately quantify gene expression changes in limited samples.

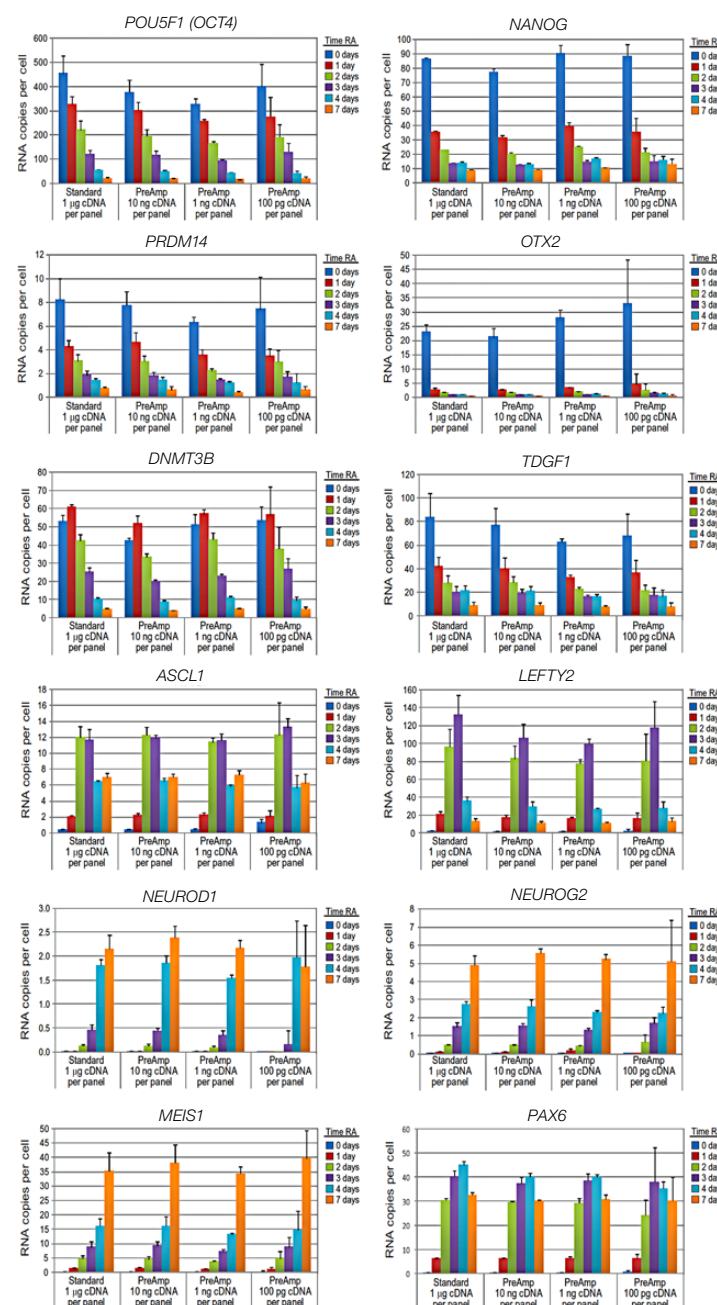


### Assessment of PreAmplification bias

NT2 cells were treated with RA for 0 to 7 days; RNA was isolated and converted to cDNA. Varied amounts of cDNA (100 pg, 1 ng and 10 ng) were pre-amplified with SsoAdvanced PreAmplification Supermix and all 100 target genes were subsequently analyzed by qPCR. In parallel, as a no-PreAmplification control, 1 µg cDNA was used to quantify the 100 gene targets directly by qPCR (10 ng/target). The experiment was performed three times independently.

The level of expression of each target gene was determined and expressed as "RNA copies per cell" after normalizing to TBP with the assumption that there are 10 TBP copies per cell. Error bars represent standard deviation of the three independent experiments. Data for target genes that show significant differentiation induced changes are shown.

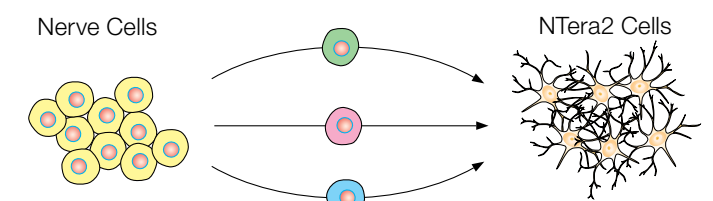
We find that, relative to standard gene expression analysis, PreAmplification does not cause statistically significant bias in the gene expression profiles of all target genes analyzed, even when 10,000-fold less cDNA is used. These findings demonstrate that SsoAdvanced PreAmplification Supermix provides accurate and reproducible quantification of both stem cell- and neural cell-specific biomarkers and does not bias gene expression profiling results.



## Single cell analysis

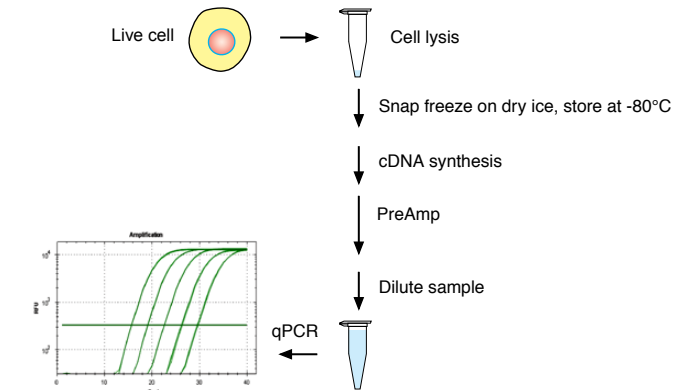
### Analysis of single cells undergoing differentiation

Most studies that address how cells change are based on population analyses. It is, however, also important to understand how individual cells behave in a changing population. For example, as a NT2 population moves from a pluripotent to a neuronal state, do individual cells change in unison, or at varying rates?



### Single cell analysis

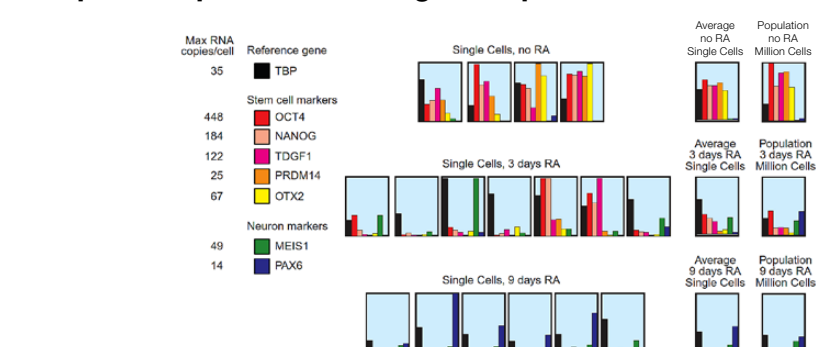
We developed a protocol that uses SsoAdvanced PreAmplification Supermix to quantify gene expression in individual NT2 cells at various time points following RA treatment. We find that cells at the 0 day time point express primarily stem cell biomarkers and cells at the 9 day time point express mainly neural cell biomarkers. Interestingly, at the 3 day time point we observe that the gene expression profiles are heterogeneous – some cells are similar to the 0 day cells while others are similar to the 9 day RA treated cells. These findings imply that NT2 cells do not undergo synchronized differentiation; instead the differentiation process is heterogeneous and cell dependent. In addition, the average gene expression profiles from the individual cells correlate with the gene expression profiles from the population analysis; this validates our single cell results.



### Gene expression in differentiating single cells

Sample	TBP	OCT4	NANOG	TDGF1	PRDM14	OTX2	MEIS1	PAX6
No cells	0	0	0	0	0	0	0	0
None Cell 1	25	131	66	69	9	9	2	0
None Cell 2	10	443	115	84	11	8	0	0
None Cell 3	23	287	104	29	25	52	0	1
None Cell 4	14	368	147	105	19	67	0	0
No cells	0	0	0	0	0	0	0	0
3d Cell 5	9	161	19	4	0	2	18	0
3d Cell 6	14	8	8	1	0	9	4	0
3d Cell 7	35	67	19	7	0	6	49	0
3d Cell 8	25	5	7	14	0	11	3	0
3d Cell 9	25	448	184	34	7	8	6	0
3d Cell 10	18	330	105	122	2	1	4	0
3d Cell 11	22	43	13	7	0	2	15	2
No Cells	0	0	0	0	0	0	0	0
9d Cell 12	6	0	1	0	0	0	5	2
9d Cell 13	18	0	1	3	0	0	4	14
9d Cell 14	10	0	0	0	0	0	4	6
9d Cell 15	6	3	1	2	0	0	2	4
9d Cell 16	10	0	5	8	0	0	5	9
9d Cell 17	19	0	5	5	0	0	9	0

### Graphical representation of gene expression in differentiating single cells



## Summary and Conclusions

We have developed a new reagent, SsoAdvanced PreAmplification Supermix, which allows for accurate quantification of up to 100 target genes from very small sample amounts including single cells. We envision that this product can benefit researchers who work with limited or rare samples, and can lead to advances in the single cell analysis field.

## References

- Andrews PW (1984). Retinoic acid induces neuronal differentiation of a cloned human embryonal carcinoma cell line in vitro. Dev Biol 103(2):285-293.
- Pleasure, SJ, and Lee, VM (1993). NTera 2 cells: a human cell line which displays characteristics expected of a human committed neuronal progenitor cell. J Neurosci Res 35: 585-602.
- Deb-Rinker, et al. (2005). Sequential DNA methylation of the Nanog and Oct-4 upstream regions in human NT2 cells during neuronal differentiation. J Biol Chem 280, 6257-60.

Use of SsoAdvanced Supermixes and PrimePCR PreAmplification Assays is covered by one or more of the following U.S. patents and corresponding patent claims outside the U.S.: 5,804,375; 5,538,848; 5,723,591; 5,876,930; 5,994,056; 6,030,787; 6,171,785; and 6,258,569. The purchase of these products includes a limited, non-transferable immunity from suit under the foregoing patent claims for using only this amount of product for the purchaser's own internal research. No right under any other patent claim and no right to perform commercial services of any kind, including without limitation reporting the results of purchaser's activities for a fee or other commercial consideration, is conveyed expressly, by implication, or by estoppel. These products are for research use only. Diagnostic uses under Roche patents require a separate license from Roche. Further information on purchasing licenses may be obtained from the Director of Licensing, Applied Biosystems, 850 Lincoln Centre Drive, Foster City, California 94404, USA.

