

SCRIPT Reverse Transcriptase

Reverse Transcriptase with increased thermal stability
Catalog-No.: PCR-505

Two-Step RT-PCR assay set-up

Reverse Transcription

Template: human HeLa cell RNA

Template dilution series: 10 ng, 1 ng, 100 pg, 10 pg

Reverse Transcription assay volume: 20 μ l

SCRIPT Reverse Transcriptase: 200 units

Primer: Oligo dT₂₀, 2.5 μ M

dNTP Mix: 500 μ M each

PCR

Assay volume: 50 μ l

Amplicon: 749 bp fragment from human GAPDH gene

Polymerase: Hot Start Pol, #PCR-203

Reverse Transcription

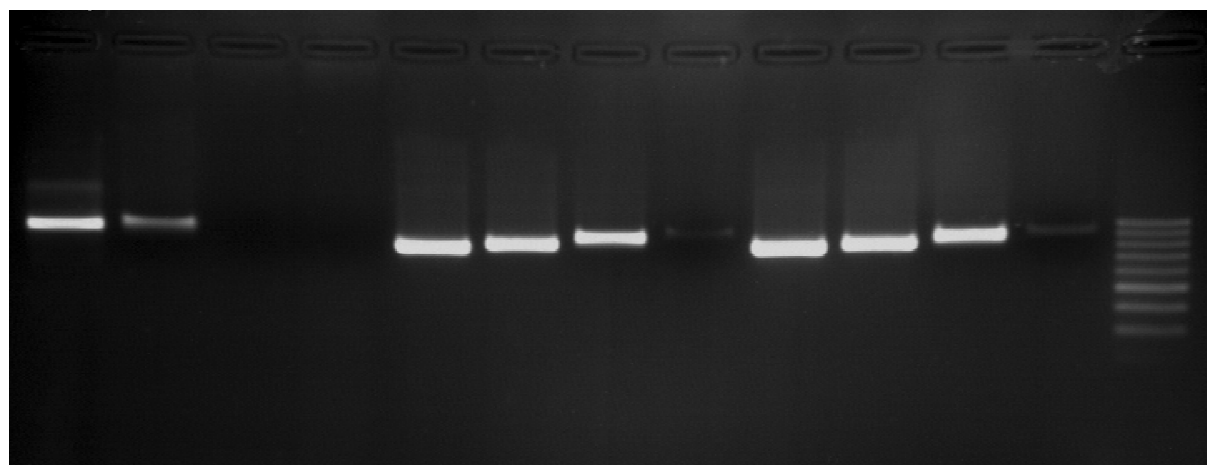
50°C	40 min	reverse transcription	1x
------	--------	-----------------------	----

PCR cycling

95°C	5 min	initial denaturation	1x
95°C	10 sec	denaturation	35x
58°C	20 sec	annealing	
72°C	1 min	elongation	
72°C	1 min	final elongation	1x

Gel Image

1,8% Agarose, 10 μ l in Gel geladen, 160 V, 30 min



10 ng	1 ng	100 pg	10 pg	10 ng	1 ng	100 pg	10 pg	10 ng	1 ng	100 pg	10 pg	100 bp Ladder
Competitor I				Competitor II				SCRIPT Reverse Transcriptase Cat.-No. PCR-505				

SCRIPT One-Step RT-PCR Kit

One-Step RT-PCR Kit for highly sensitive and specific amplification
 Catalog-No.: PCR-509

One-Step RT-PCR assay set-up

Template: human HeLa cell RNA

Template dilution series: 10 ng, 1 ng, 100 pg, 10 pg

Reverse Transcription assay volume: 50 μ l

Primer: 400 nM each

Amplicon: 749 bp fragment from human GAPDH gene

RT-PCR cycling

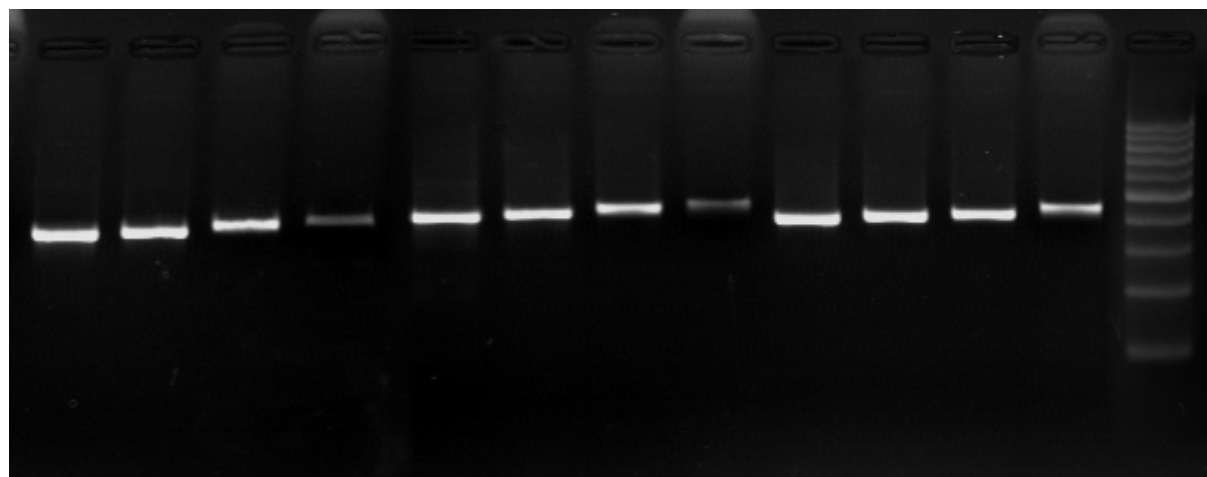
50°C	40 min	reverse transcription	1x
95°C	2 min	initial denaturation	1x
95°C	10 sec	denaturation	35x
58°C	20 sec	annealing	
72°C	1 min	elongation	
72°C	1 min	final elongation	1x

Gel Image

1,8% Agarose

10 μ l in Gel geladen

160 V, 30 min



10 ng	1 ng	100 pg	10 pg	10 ng	1 ng	100 pg	10 pg	10 ng	1 ng	100 pg	10 pg	200 bp Ladder
Competitor I				Competitor II				SCRIPT One-Step RT-PCR Kit Cat.-No. PCR-509				

SCRIPT High Fidelity One-Step RT-PCR Kit

One-Step RT-PCR Kit for highly precise and fast amplification
Catalog-No.: PCR-510

One-Step RT-PCR assay set-up

Template: human HeLa cell RNA

Template dilution series: 10 ng, 1 ng, 100 pg, 10 pg

Reverse Transcription assay volume: 50 μ l

Primer: 400 nM each

Amplicon: 749 bp fragment from human GAPDH gene

RT-PCR cycling

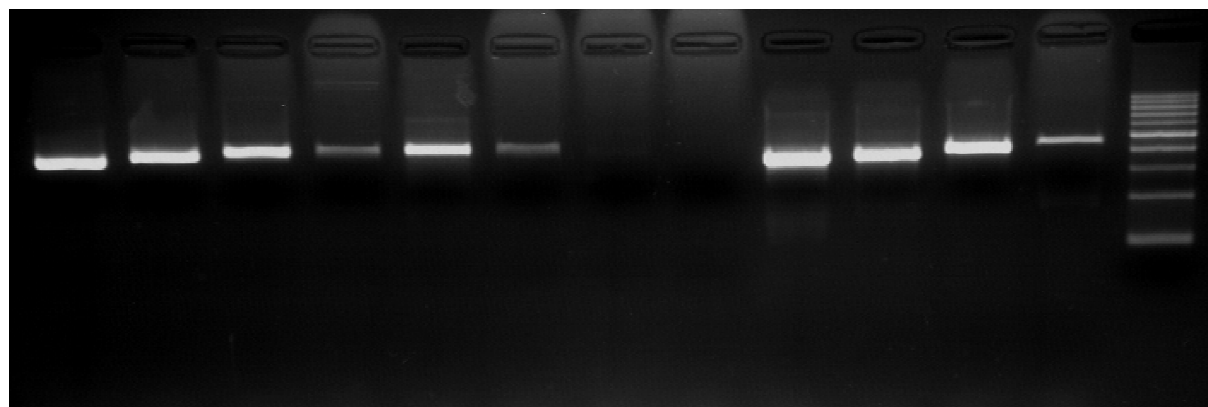
50°C	40 min	reverse transcription	1x
95°C	2 min	initial denaturation	1x
95°C	10 sec	denaturation	35x
58°C	20 sec	annealing	
72°C	30 sec	elongation	
72°C	1 min	final elongation	1x

Gel Image

1,8% Agarose

10 μ l in Gel geladen

160 V, 30 min



10 ng	1 ng	100 pg	10 pg	10 ng	1 ng	100 pg	10 pg	10 ng	1 ng	100 pg	10 pg	200 bp Ladder
Competitor I				Competitor II				SCRIPT High Fidelity One-Step RT-PCR Kit Cat.-No. PCR-510				