

THE PCR SONG

Words and Music by
Tyler Kay

Arranged by
Mac Huff

BIO - RAD LABORATORIES

BIO-RAD

The PCR Song

There was a time when to amplify DNA,
You had to grow tons and tons of tiny cells.

Then along came a guy named Dr. Kary Mullis,
Said you could amplify *in vitro* just as well.

Just mix your template with a buffer & some primers ~
Nucleotides and polymerases, too.

Denaturing, annealing and extending.
Well it's amazing what heating and cooling and heating will do.

PCR, when you need to detect mutations.

PCR, when you need to recombine.

PCR, when you need to find out who the daddy is.

PCR, when you need to solve a crime.

(repeat chorus)



The PCR Song

Words and Music by
Tyler Kay
Arranged by Mac Huff

Strong Anthem Rock

Lead

Bk-up

S/A

T/B

There was a time when to am-pl-fy D. N. A. You

solo 1

Strong Anthem Rock

Piano

80's rhodes dbl.

G C G/B Ami7 G C D G D/F# Emi

Strong Anthem Rock

Synth.

Strong Anthem Rock

Guitar

C D G D/F# Emi

Strong Anthem Rock

Bass

C D G D/F# Emi

Strong Anthem Rock

Drums

pre roll - mallets 2 3 4 (Rim Knocks)

Lead

Bk-up

S/A

T/B

Piano

Synth.

Guitar

Bass

Drums

had to grow tons and tons of ti - ny cells. Then a - long came a guy named Doctor Ka - ry Mul - lis 'said you can

solo 2

solo 3

Ami G D/F# D D7/C Bmi7 Ami7 G D/F# D

Ami G D/F# D D7/C Bmi7 Ami7 G D/F# D

5 6 7 8

Detailed description: This is a musical score for a song, likely a pop or rock track. It consists of nine staves. The top two staves are for Lead and Bk-up, both of which are currently blank. The third staff is for S/A (Soprano/Alto) and contains the lyrics: "had to grow tons and tons of ti - ny cells. Then a - long came a guy named Doctor Ka - ry Mul - lis 'said you can". Above the lyrics, there are musical notations including eighth and sixteenth notes, rests, and two solo markings labeled "solo 2" and "solo 3". The fourth staff is for T/B (Tenor/Bass), which is also blank. The fifth staff is for Piano, showing a bass line with chords and a treble line with chords and some melodic movement. Above the piano staff, a series of chords are listed: Ami, G, D/F#, D, D7/C, Bmi7, Ami7, G, D/F#, D. The sixth staff is for Synth., which is blank. The seventh staff is for Guitar, showing a rhythmic pattern with slanted lines and some chordal movement. Above the guitar staff, the same series of chords as above the piano staff are listed: Ami, G, D/F#, D, D7/C, Bmi7, Ami7, G, D/F#, D. The eighth staff is for Bass, showing a bass line with notes and rests. Above the bass staff, the same series of chords as above the piano staff is listed: Ami, G, D/F#, D, D7/C, Bmi7, Ami7, G, D/F#, D. The ninth staff is for Drums, showing a rhythmic pattern with slanted lines. At the bottom of the drum staff, the numbers 5, 6, 7, and 8 are written, corresponding to the measures.



Lead

Bk-up

S/A
am pli fy in vi - tro just as well. Just mix your tem - plate with a buffer and some pri - mers.

T/B

Piano
Ami7 G/B C D G D/F# Emi

Synth.

Guitar
Ami7 G/B C D G D/F# Emi

Bass
Ami7 G/B C D G D/F# Emi

Drums
Fill

9 10 11 12

Lead

Bk-up

S/A

T/B

Piano

Synth.

Guitar

Bass

Drums

oo

solo 5

solo 6

Nu-cle - o - tides and po - ly - mer - a - ses too. De - nat - ur - ing an - neal - ing and ex - tend ing Well it's a -

Ami G D/F# D D7/C Bmi7 Ami7 G D/F# D

Ami G D/F# D D7/C Bmi7 Ami7 G D/F# D

13 14 15 16

19

Lead

P C R when you need to de tect mu ta tions, P C

Bk-up

oo Ah de - tect mu - ta - tions

S/A

mazing what heating and cooling and heating will do. P C R when you need to de tect mu ta tions, C

T/B

Piano

Ami7 G/B C D C/D 19 G D/F# Emi

Synth.

Guitar

Ami7 G/B C D C/D 19 G D/F# Emi

add pwr. chrd

Bass

Ami7 G/B C D C/D 19 G D/F# Emi

Drums

Fill _____ 19 to big sn.

17 18 19 20

Lead
R when you need to re - com-bine. P C R when you need to find out who the

Bk-up
Ah re - com-bine. Ah

S/A
R when you need to re - com-bine. P C R when you need to find out who the

T/B

Piano
Ami G D/F# D D7/C Bmi7 C

Synth.

Guitar
Ami G D/F# D D7/C Bmi7 C

Bass
Ami G D/F# D D7/C Bmi7 C

Drums
Fill

21 22 23

Lead
dad - dy is, P C R when you need to solve a crime. P C

Bk-up
Who's your dad dy. Ah solve a crime.

S/A
dad - dy is, P C R when you need to solve a crime. P C

T/B

Piano
G/D D Ami7 G/B D A D/F#

Synth.

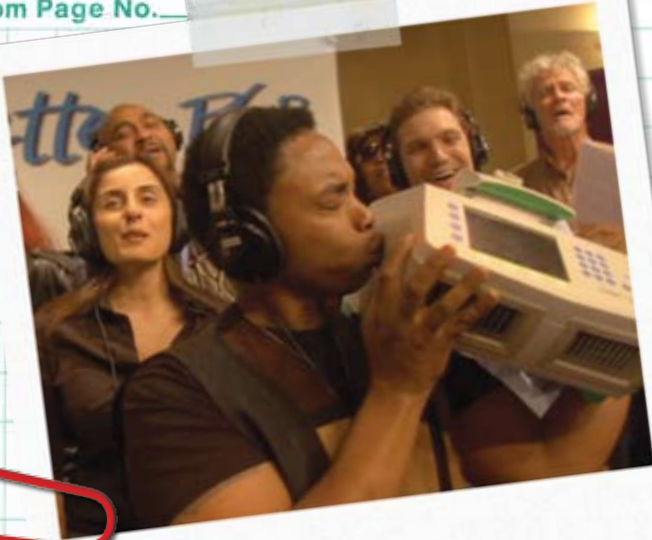
Guitar
G/D D Ami7 G/B D A D/F#

Bass
G/D D Ami7 G/B D A D/F#

Drums
24 25 26

TITLE _____

From Page No. _____



To Page N

Witnessed & Understood by me,

Date

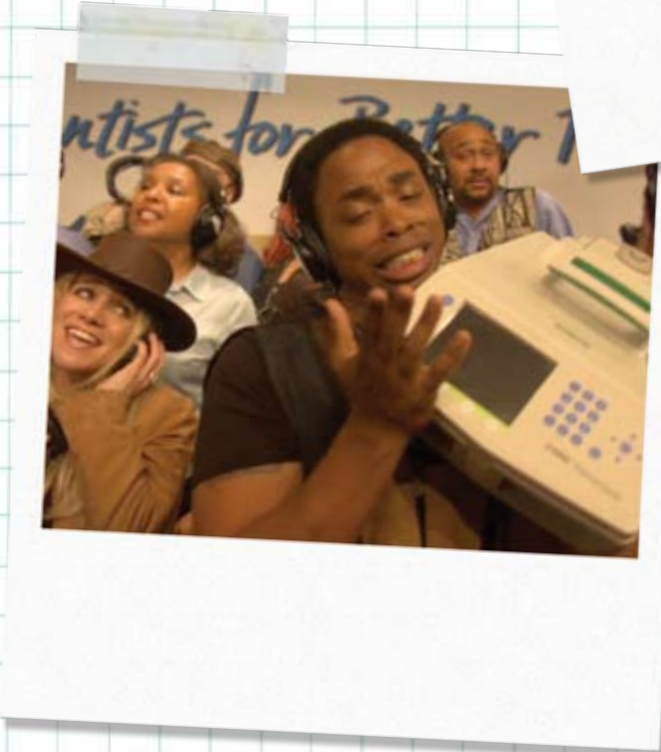
Invented by

Date

Recorded by

TITLE _____

From Page _____



Practice of polymerase chain reaction (PCR) may require a licence.

To Page N

Witnessed & Understood by me,

Date

Invented by

Date

Recorded by



**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Web site www.bio-rad.com **USA** 800 4BIORAD **Australia** 61 02 9914 2800 **Austria** 01 877 89 01 **Belgium** 09 385 55 11 **Brazil** 55 21 3237 9400
Canada 905 364 3435 **China** 86 21 6426 0808 **Czech Republic** 420 241 430 532 **Denmark** 44 52 10 00 **Finland** 09 804 22 00 **France** 01 47 95 69 65
Germany 089 318 84 0 **Greece** 30 210 777 4396 **Hong Kong** 852 2789 3300 **Hungary** 36 1 455 8800 **India** 91 124 4029300 **Israel** 03 963 6050
Italy 39 02 216091 **Japan** 03 6361 7000 **Korea** 82 2 3473 4460 **Mexico** 52 555 488 7670 **The Netherlands** 0318 540666 **New Zealand** 0508 805 500
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** 061 717 95 55 **Taiwan** 886 2 2578 7189 **United Kingdom** 020 8328 2000