

GENE TRANSFER

Gene Pulser MXcell™ Electroporation System

Preset Protocol Quick Guide

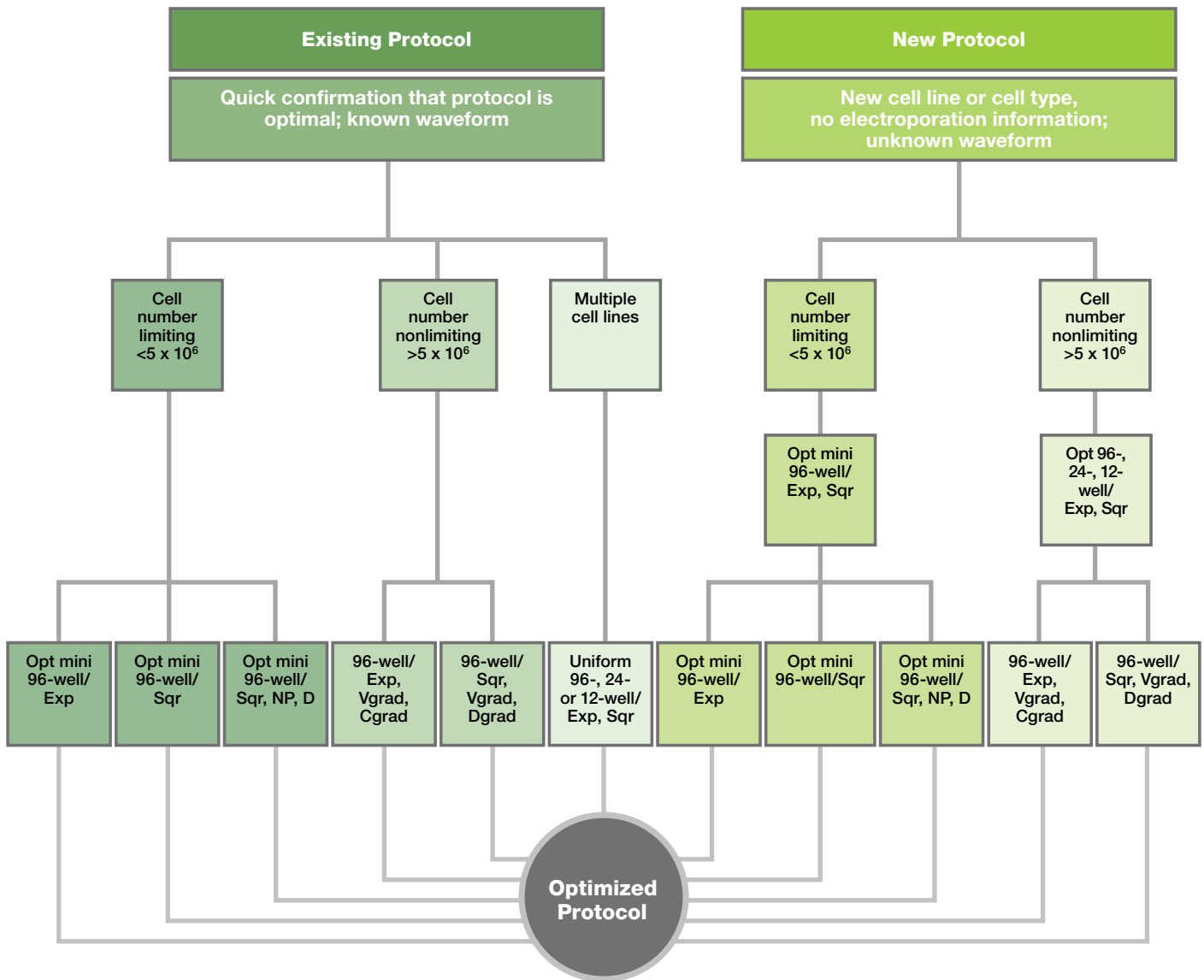


Preset Protocol*	Plate	Use
Opt mini 96-well/Sqr, Exp	96	To rapidly determine optimal waveform and conditions
Opt mini 96-well/Sqr	96	To rapidly determine optimal conditions for square-wave protocols
Opt mini 96-well/Exp	96	To rapidly determine optimal conditions for exponential protocols
Opt 96-well/Sqr, NP, D	96	To enhance cell viability and improve efficiency after optimal square-wave conditions have been determined
96-well/Exp	96	For initial protocol setup with exponential waveform conditions
24-well/Exp	24	For initial protocol setup with exponential waveform conditions
12-well/Exp	12	For initial protocol setup with exponential waveform conditions
96-well/Sqr	96	For initial protocol setup with square-wave conditions
24-well/Sqr	24	For initial protocol setup with square-wave conditions
12-well/Sqr	12	For initial protocol setup with square-wave conditions
96-well/Exp, Vgrad, Cgrad	96	To fine-tune conditions with new or existing cell lines when exponential waveforms are normally applied; includes replicates
96-well/Sqr, Vgrad, Dgrad	96	To fine-tune conditions with new or existing cell lines when square waveforms are normally applied; includes replicates
Opt 96-well/Exp, Sqr	96	For cell lines with no protocol reference; includes a range of average starting conditions
Opt 24-well/Exp, Sqr	24	For cell lines with no protocol reference; includes a range of average starting conditions
Opt 12-well/Exp, Sqr	12	For cell lines with no protocol reference; includes a range of average starting conditions
Uniform 96-well/Exp, Sqr	96	To compare different cell lines or electroporation of different molecules within or between cell lines
Uniform 24-well/Exp, Sqr	24	To compare different cell lines or electroporation of different molecules within or between cell lines
Uniform 12-well/Exp, Sqr	12	To compare different cell lines or electroporation of different molecules within or between cell lines
Mixed 96-well/Exp, Sqr	96	For mixing different waveforms by alternating rows of exponential (250 V/350 μ F) and square waveforms (250 V/20 ms)
Mixed 24-well/Exp, Sqr	24	For mixing different waveforms by alternating rows of exponential (250 V/350 μ F) and square waveforms (250 V/20 ms)
Mixed 12-well/Exp, Sqr	12	For mixing different waveforms by alternating rows of exponential (250 V/350 μ F) and square waveforms (250 V/20 ms)

* Protocol nomenclature — First word describes protocol purpose (Opt, optimal), followed by electroporation plate type (for example, mini 96-well), followed by slash and waveform type (Exp, exponential; Sqr, square-wave), followed by specific parameters, if applicable (V, voltage; R, resistance; C, capacitance; D, pulse duration; NP, number of pulses; Grad, gradient). Note: General protocols do not include protocol purpose.



Preset Protocol Decision Tree



**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 905 712 2771 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 5811 6270 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