

ddPCR™ *SMN1* Copy Number Determination Kit



The ddPCR *SMN1* Copy Number Determination Kit is designed to determine the copy number of the *SMN1* gene only. The kit contains a 20x ddPCR *SMN1* Copy Number Determination Assay, positive controls, and ddPCR Supermix for Probes (No dUTP).

Combined with Bio-Rad's QX100™, QX200™, or QX200™ AutoDG™ Droplet Digital PCR (ddPCR) System, the ddPCR *SMN1* Kit allows you to:

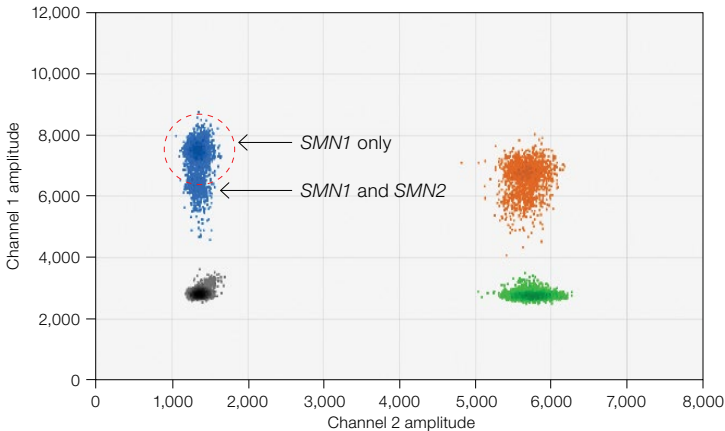
- **Generate accurate and reproducible copy number calls for *SMN1*** — provides superior performance
- **Perform high-throughput screening** — allows processing of several hundred samples in a single day
- **Benefit from economical cost per sample**

Visit bio-rad.com/web/SMN1CNkit for more information.

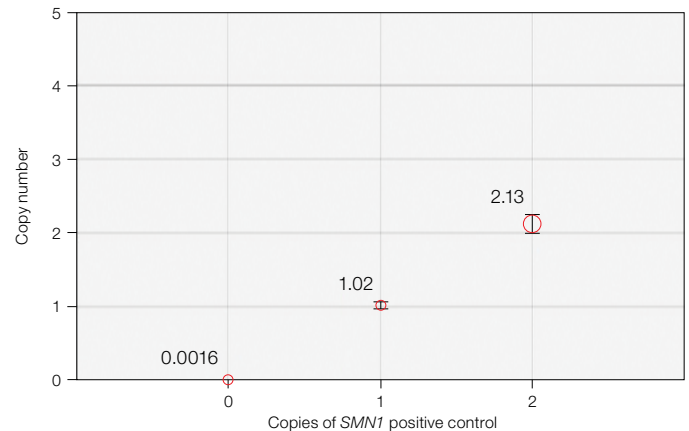
BIO-RAD

Near-Integer Copy Number Calls for *SMN1*

Spinal muscular atrophy (SMA) is a group of inherited diseases that cause muscles to lose function. The progressive muscle deterioration causes weakness and eventually leads to death. SMA affects the motor neurons, especially those in the spinal cord and brain stem. The gene affected is known as the survival motor neuron gene (*SMN1*) and is normally responsible for creating a protein necessary for motor neurons to function correctly. The ddPCR *SMN1* Copy Number Determination Kit is designed to detect the copy number of the *SMN1* gene only. For *SMN2* detection, use the ddPCR *SMN2* Copy Number Determination Kit (catalog #186-3503).



Detection of *SMN1* copy number. A 2-D fluorescence amplitude plot shows the wells of a sample duplexed with the *SMN1* and reference assays. The black cluster on the plot represents the negative droplets, the green cluster represents the droplets that are positive for the reference only, the blue cluster within the red circle represents the droplets that are present for *SMN1* only, and the orange cluster represents the droplets that are positive for both *SMN1* and the reference target.



Accurate and reproducible copy number calls for *SMN1*. Copy number plot shows positive controls, ranging from zero to two copies, detected using the ddPCR *SMN1* Copy Number Determination Kit. All error bars are generated by QuantaSoft™ Software and represent a 95% confidence interval.

Ordering Information

Catalog #	Description
186-3500	ddPCR <i>SMN1</i> Copy Number Determination Kit, 200 x 20 µl reactions, includes 20x ddPCR <i>SMN1</i> Copy Number Determination Assay (1 x 200 µl vial), positive controls, ddPCR Supermix for Probes (No dUTP) (2 x 1 ml vials)

The QX100 or QX200 Droplet Digital PCR System and/or its use is covered by claims of U.S. patents, and/or pending U.S. and non-U.S. patent applications owned by or under license to Bio-Rad Laboratories, Inc. Purchase of the product includes a limited, non-transferable right under such intellectual property for use of the product for internal research purposes only. No rights are granted for diagnostic uses. No rights are granted for use of the product for commercial applications of any kind, including but not limited to manufacturing, quality control, or commercial services, such as contract services or fee for services. Information concerning a license for such uses can be obtained from Bio-Rad Laboratories. It is the responsibility of the purchaser/end user to acquire any additional intellectual property rights that may be required.



**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Web site www.bio-rad.com USA 800 424 6723 Australia 61 2 9914 2800 Austria 43 1 877 89 01 Belgium 03 710 53 00 Brazil 55 11 3065 7550
Canada 905 364 3435 China 86 21 6169 8500 Czech Republic 420 241 430 532 Denmark 44 52 10 00 Finland 09 804 22 00
France 01 47 95 69 65 Germany 49 89 31 884 0 Greece 30 210 9532 220 Hong Kong 852 2789 3300 Hungary 36 1 459 6100 India 91 124 4029300
Israel 03 963 6050 Italy 39 02 216091 Japan 81 3 6361 7000 Korea 82 2 3473 4460 Mexico 52 555 488 7670 The Netherlands 0318 540666
New Zealand 64 9 415 2280 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 (0) 861 246 723 Spain 34 91 590 5200 Sweden 08 555 12700 Switzerland 026674 55 05
Taiwan 886 2 2578 7189 Thailand 1800 88 22 88 United Kingdom 020 8328 2000