



## Droplet Digital™ PCR (ddPCR™) Supermixes



Bio-Rad's ddPCR supermix for probes and ddPCR supermix for probes (no dUTP) are ready-to-use 2x universal mixes that have been optimized to deliver maximum PCR efficiency, specificity, and sensitivity in Droplet Digital PCR (ddPCR). Both supermixes contain all the components, except primers, probe(s), and template, required for hydrolysis probe-based ddPCR.

Combined with Bio-Rad's Droplet Digital PCR system, both supermixes allow you to:

- Amplify and detect targets using commercially available probe-based assays
- Amplify from a wide range of input samples, including genomic DNA, cDNA, and plasmid
- Obtain precise and sensitive absolute measurements for a variety of applications, such as copy number analysis and mutation detection

For more information, visit [www.bio-rad.com/web/ddPCRSupermixProbes](http://www.bio-rad.com/web/ddPCRSupermixProbes).

**BIO-RAD**

## Optimized for Probe-Based ddPCR

Both supermixes have been optimized to deliver maximum PCR efficiency, specificity, and sensitivity for probe-based ddPCR. ddPCR supermix for probes is suitable for use with uracil N-glycosylase (UNG) decontamination protocols, which prevent the reamplification of carryover PCR products between experiments.

ddPCR supermix for probes (no dUTP) is highly recommended for use with all of Bio-Rad's pre-designed and fully wet-lab validated PrimePCR™ ddPCR assays and is also suitable for applications such as library quantification and sample preparation for next-generation sequencing. The table highlights the recommended applications and considerations for each supermix.



Applications and Considerations	ddPCR Supermix for Probes	ddPCR Supermix for Probes (no dUTP)
Suitable for UNG decontamination protocols	√	—
Compatible with validated PrimePCR ddPCR mutation detection assays	•	√
Compatible with validated PrimePCR ddPCR copy number assays	•	√
ddPCR library quantification kit for Illumina TruSeq	—	√
ddPCR library quantification kit for Ion Torrent	—	√
Amplification in droplets for downstream sequencing	—	√

√ Recommended supermix

• Compatible supermix

## Ordering Information

Catalog #	Description
186-3026	<b>ddPCR Supermix for Probes</b> , 2 ml (2 x 1 ml), 200 x 20 µl reactions, 2x supermix
186-3010	<b>ddPCR Supermix for Probes</b> , 5 ml (5 x 1 ml), 500 x 20 µl reactions, 2x supermix
186-3027	<b>ddPCR Supermix for Probes</b> , 25 ml (5 x 5 ml), 2,500 x 20 µl reactions, 2x supermix
186-3028	<b>ddPCR Supermix for Probes</b> , 50 ml (10 x 5 ml), 5,000 x 20 µl reactions, 2x supermix
186-3023	<b>ddPCR Supermix for Probes (no dUTP)</b> , 2 ml (2 x 1 ml), 200 x 20 µl reactions, 2x supermix
186-3024	<b>ddPCR Supermix for Probes (no dUTP)</b> , 5 ml (5 x 1 ml), 500 x 20 µl reactions, 2x supermix
186-3025	<b>ddPCR Supermix for Probes (no dUTP)</b> , 25 ml (5 x 5 ml), 2,500 x 20 µl reactions, 2x supermix

Illumina and TruSeq are trademarks of Illumina, Inc. Illumina is not affiliated with Bio-Rad Laboratories, Inc. Ion Torrent is a trademark of Life Technologies Corporation.



**Bio-Rad  
Laboratories, Inc.**

Life Science  
Group

**Web site** [www.bio-rad.com](http://www.bio-rad.com) **USA** 800 424 6723 **Australia** 61 2 9914 2800 **Austria** 01 877 89 01 **Belgium** 09 385 55 11 **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** 089 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** 27 861 246 723 **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