



**Bio-Rad
Laboratories, Inc.**

Digital Biology Center
5731 West Las Positas Blvd.
Pleasanton, California 94588
Phone: 925-474-8600
Fax: 925-474-8644

August 1, 2017

Subject: ddPCR diagnostics Product Brand Name change

Dear Valued Customer,

Since the initial launch of Bio-Rad Droplet Digital PCR platform, Bio-Rad has been using QX200™ Droplet Digital PCR to identify its systems brand.

As Bio-Rad is expanding the use of its ddPCR platform to in vitro diagnostics (IVD), we have decided to identify the diagnostics product line with the new brand name: **QXDx™**. Hence, we will be updating any current and future diagnostics digital PCR product with the prefix **QXDx™**.

The product name change does not affect the form, fit, or function, or claims of the specific products.

Starting June 15th, 2017, documentation will be updated in a phased process to reflect the **QXDx™** brand name in all diagnostics digital PCR products.

During the transition and until we complete registration of the new name, we will use both the current registered name and the new proprietary name to refer to the instruments and products as shown below:

Instrumentation Systems:

Catalog #	Registered Name	New Proprietary/Common Name
17002229	QX200™ AutoDG™ Droplet Digital PCR Dx System	QXDx™ AutoDG™ ddPCR™ System
12001045	QX200™ Droplet Reader, IVD	QXDx™ Droplet Reader
12001630	QX200™ Automated Droplet Generator, IVD	QXDx™ Automated Droplet Generator
17000034	QX200™ Droplet Digital PCR Dx System	QXDx™ ddPCR™ System
12001045	QX200™ Droplet Reader, IVD	QXDx™ Droplet Reader
12001049	QX200™ Droplet Generator, IVD	QXDx™ Droplet Generator

Reagents & Consumables:

Catalog #	Description	New Name
17001378	ddPCR™ Dx Universal Kit for AutoDG™	QXDx™ Universal Kit for AutoDG™ ddPCR™ System

12001922	ddPCR™ Dx AutoDG Consumable Pack	QXDx™ AutoDG™ Consumable Pack
12003031	ddPCR™ Dx AutoDG Supermix Pack	QXDx™ AutoDG™ Supermix Pack
12002526	ddPCR™ Dx Droplet Reader Oil Pack	QXDx™ Droplet Reader Oil Pack
17001379	ddPCR™ Dx Universal Kit for QX200™ ddPCR System	QXDx™ Universal Kit for ddPCR™ System
12001921	ddPCR™ Dx Consumable Pack	QXDx™ Consumable Pack
12002544	ddPCR™ Dx Supermix Pack	QXDx™ Supermix Pack
12002526	ddPCR™ Dx Droplet Reader Oil Pack	QXDx™ Droplet Reader Oil Pack

Sincerely,

Marketing Manager,
Bio-Rad, Digital Biology Center