

## Important Note

Roche discontinuation of LightCycler TaqMan Master (Master Mix PCR 5x) (Roche, cat. no. 04535286001) and LightCycler FastStart DNA MasterPLUS HybProbe® (Master Mix 5x) (Roche, cat. No. 03515567001)

Dear valued customer,

You may have been informed by Roche about the discontinuation effective on 31<sup>st</sup> December 2024 of their LightCycler TaqMan Master (Master Mix PCR 5x) (Roche, cat. no. 04535286001) and LightCycler FastStart DNA MasterPLUS HybProbe® (Master Mix 5x) (Roche, cat. No. 03515567001).

This discontinuation solely affects the *ipsogen*® products (listed in below table), used with Roche real-time qPCR capillary instruments LightCycler 1.2, LightCycler 1.5, and LightCycler 2.0 that are no longer supported, or will no longer be supported by Roche after 31<sup>st</sup> December 2024.

If you are affected by the discontinuation, you can still use the listed *ipsogen* products with alternative non-capillary real-time qPCR instruments and alternative reagents listed in the “Materials Required but Not Provided” section of the respective handbook, following their related instructions for use.

| Affected <i>ipsogen</i> product Cat. No. | Affected <i>ipsogen</i> product Name       |
|------------------------------------------|--------------------------------------------|
| 670023                                   | <i>ipsogen</i> BCR-ABL1 mbcr Kit (24) CE   |
| 670123                                   | <i>ipsogen</i> BCR-ABL1 Mbcr Kit (24) CE   |
| 670125                                   | <i>ipsogen</i> BCR-ABL1 Mbcr Kit CE (52)   |
| 673522                                   | <i>ipsogen</i> JAK2 MutaQuant Kit (12) CE  |
| 673523                                   | <i>ipsogen</i> JAK2 MutaQuant Kit (24) CE  |
| 673022                                   | <i>ipsogen</i> JAK2 MutaScreen Kit (10) CE |
| 673023                                   | <i>ipsogen</i> JAK2 MutaScreen Kit (24) CE |

|        |                                                |
|--------|------------------------------------------------|
| 673123 | <i>ipsogen</i> JAK2 MutaScreen RS Kit (19) CE  |
| 673823 | <i>ipsogen</i> JAK2 MutaSearch Kit (24) CE     |
| 672123 | <i>ipsogen</i> PML-RARA bcr1 Kit (24) CE       |
| 676923 | <i>ipsogen</i> WT1 ProfileQuant Kit (24) CE    |
| 670013 | <i>ipsogen</i> BCR-ABL1 mbcr Kit (24)          |
| 670113 | <i>ipsogen</i> BCR-ABL1 Mbcr Kit (24)          |
| 670115 | <i>ipsogen</i> BCR-ABL1 Mbcr Kit (52)          |
| 672113 | <i>ipsogen</i> PML-RARA bcr1 Kit (24)          |
| 672213 | <i>ipsogen</i> PML-RARA bcr2 Kit (24)          |
| 672313 | <i>ipsogen</i> PML-RARA bcr3 Kit (24)          |
| 675013 | <i>ipsogen</i> RUNX1-RUNX1T1 Kit (24)          |
| 675113 | <i>ipsogen</i> ETV6-RUNX1 Kit (24)             |
| 676013 | <i>ipsogen</i> CBFβ-MYH11 A Kit (24)           |
| 673512 | <i>ipsogen</i> JAK2 MutaQuant Kit (12)         |
| 673513 | <i>ipsogen</i> JAK2 MutaQuant Kit (24)         |
| 673013 | <i>ipsogen</i> JAK2 MutaScreen Kit (24)        |
| 673012 | <i>ipsogen</i> JAK2 MutaScreen Kit (10)        |
| 673113 | <i>ipsogen</i> JAK2 MutaScreen RS Kit (19)     |
| 673813 | <i>ipsogen</i> JAK2 MutaSearch Kit (24)        |
| 676413 | <i>ipsogen</i> MPL W515L/K MutaScreen Kit (24) |
| 676913 | <i>ipsogen</i> WT1 ProfileQuant Kit (24)       |
| 676613 | <i>ipsogen</i> BAALC ProfileQuant Kit (24)     |
| 677513 | <i>ipsogen</i> NPM1 mut A MutaQuant Kit (24)   |
| 677613 | <i>ipsogen</i> NPM1 mut B&D MutaQuant Kit (24) |

We very much value our relationship with you and your laboratory, and we apologize for any inconvenience this Roche discontinuation may cause. We hope that you will find the proposed alternative useful.

With kind regards,

QIAGEN

[www.qiagen.com](http://www.qiagen.com)