

June 2021

RE: Discontinuation of 5 ml tube for PCR mix and new catalog number announcement

Dear Valued Customer,

We would like to inform you about the discontinuation of the product listed below. This change is effective immediately. This product will no longer be manufactured and will be available until June 2021 or until the current inventory is depleted.

Catalog #	Discontinued Product	Alternative Catalog #	Product Name
3594901	5 ml Tube for PCR Mix (50 tubes/unit)	12016673	5 ml Tube for PCR Mix (25 tubes/unit) (available for order starting June 18, 2021)

The new 5 ml tube for PCR mix has been validated to work on the iQ-Check Prep System. The main changes to the new item are listed as follows:

- The current 5 ml tube for PCR mix does not include a cap on the tube. The new 5 ml tube for PCR mix will come with a white cap on the tube.
- The current 5 ml tube for PCR mix are packaged in 50 tubes per bag. The new 5 ml tube for PCR mix will be packaged in 25 tubes per bag.

We apologize for any inconvenience this may cause. Please do not hesitate to contact your Bio-Rad representative with any questions.

Best Regards,

Jean-Philippe Tourniaire
International Product Manager
Bio-Rad Laboratories, Food Science Division