



Date: February 7th 2022
From: Randi Muster, Global Commercial Product Manager
Re: Allegra X-5 Obsolescence

Dear Valued Customer,

Beckman Coulter Life Sciences is dedicated to providing you with the highest quality support and communicating information regarding our products. As a part of our continuous improvement procedures, we periodically evaluate our product portfolio to optimize our offerings. Based on this evaluation, we are obsoleting our Allegra X-5 benchtop centrifuge and will replace it with the **Allegra V-15R**.

The Allegra X-5 products listed below will no longer be available **January 31, 2022**.

We will continue to fully service and support the Allegra X-5 centrifuge for a period of five (5) years after the obsolescence date (i.e., through February 28, 2027). During this time, routine preventive maintenance, repairs, service agreements, and all other support activities will be available.

Obsoleted Allegra X-5 Models Effective January 31, 2021:

Part No.	Description
B30585	Allegra X-5 IVD 120V 60 Hz
B30586	Allegra X-5 IVD 220V-240V 50/60 Hz
B30587	Allegra X-5 IVD 200V 50/60 Hz
B30588	Allegra X-5 IVD 220V-240V w/CCC plug
B30589	Allegra X-5 Clinical Sample Prep Package, 120 V 60 Hz, IVD
B30590	Allegra X-5 Clinical Sample Prep Package, 220-240 V 50-60 Hz, IVD
B30591	Allegra X-5 Clinical Sample Prep Package, 200 V 50-60 Hz, IVD
B30592	Allegra X-5 Clinical Sample Prep Package, 220-240 V w/CCC plug, IVD

Allegra V-15R Benchtop Centrifuge

The Allegra V-15R Benchtop Centrifuge is a versatile workhorse compatible with 10 available rotors, which enables numerous configurations and allow you to spin up to 750 mL bottles, various tubes and microplates.

High Level Features:

- Compact footprint, which allows this centrifuge to easily fit on a lab bench
- Enhanced safety features such as automatic rotor recognition and imbalance sensors
- Easy-to-use interface with illuminated button to quickly display run status

Comparison Chart:

Feature	Allegra V-15R	Allegra X-5
Max RPM	13,500	4,700
Max RCF	20,412 x g	4,470 x g
Max Capacity	4 x 750 mL	4 x 100 x 10 mL
50 mL Conical Tubes	28	--
15 mL Conical Tubes	60	--
10 mL RIA	100	100
5/7 mL RIA	144	140
No. of Rotors	10	1

If you'd like more information regarding replacement options, please contact us at www.beckman.com/contact-us

We look forward to continuing to provide you with all the centrifuges, rotors and labware that you need for your research.

Sincerely,

Randi Muster

Randi Muster

Global Commercial Product Manager, *Beckman Coulter Life Sciences*
rmuster@beckman.com