



Date: November 30, 2021
From: Randi Muster, Global Commercial Product Manager
Re: Avanti J-XPI/XP 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 product offerings. Based on this evaluation, we are obsoleting our Avanti J-XPI/XP high performance centrifuges as there are new models available.

The Avanti J-I/XPI/XP products listed below will no longer be available for purchase as of **January 31, 2022**.

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

Obsoleted Avanti J-I/XPI/XP Models Effective January 31, 2022:

- ***Avanti J-30I***
- ***Avanti J-26XP***
- ***Avanti J-26XPI***

Avanti JXN High Performance Centrifuges

The Avanti JXN Series centrifuges are designed for high throughput and productivity and are ideal for shared labs and GMP environments requiring enhanced biocontainment and advanced levels of networking and data management.

High Level Features:

- GMP Supporting software through large, intuitive user interface
- Monitoring and controlled flexibility with MobileFuge
- Enhanced safety features such as automatic rotor recognition and imbalance sensors

Avanti Model Comparison Chart:

Feature	JXN-30	JXN-26	J-30i	J-26XPI	J-26XP
Max RPM	30,000	26,000	30,000	26,000	26,000
Max RCF	110,500 x g	81,770 x g	110,500 x g	81,800 x g	81,000 x g
Max Capacity	4 L	6 L	4 L	4 L	4 L
GMP Supporting	Yes	Yes	--	--	--
Interface	15" Color Touchscreen LCD	15" Color Touchscreen LCD	Backlit Display w/ Keypad	Backlit Display w/ Keypad	Backlit Display w/ Knobs

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

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