

Sartorius Lab Instruments GmbH & Co. KG, 37070 Göttingen

Changenotification_ab@sartorius.com

October 26, 2023

Expected implementation date	Customer action required by Sartorius	Type of action
January 2024	No	N/A

Notification Letter

ECR 41686: Discontinuation of Picus® and Picus®Nxt Pipettes

Dear Valued Customer,

This change notification is to inform you of the discontinuation of Picus® and Picus® Nxt electronic pipettes.

Change Description

As part of our life-cycle management process, we are streamlining our offerings. The recently released Picus® 2 electronic pipette will serve as a replacement for both, the Picus® and Picus® Nxt electronic pipettes.

Impacted Products

The list of impacted products can be found in Appendix 1 at the end of this letter.

Change Implementation

This change is controlled internally in accordance with Sartorius change management procedures.

The product manufacturing will be stopped in January 2024. To ensure availability, we recommend placing your final orders no later than December 31, 2023. Please contact Sartorius customer service for larger final orders. Sartorius distribution centers will serve customer orders as long as stocks last after discontinuation.

Picus® and Picus® Nxt electronic pipettes are replaced by Picus® 2 electronic pipettes. For more information about alternative products, please refer to Appendix 1, consult your Sartorius Sales representative, or visit <https://www.sartorius.com/en/products/pipetting/electronic-pipettes/picus-2>

Sartorius Lab Instruments GmbH & Co. KG

Otto-Brenner-Straße 20
37079 Göttingen, Germany
Phone +49 551 308 0
Fax +49 551 308 3289
www.sartorius.com

Registered Office:
Göttingen
Local Court of Registration:
Amtsgericht Göttingen
HRA 201253

The company is a limited
commercial partnership

Individually liable shareholder

Sartorius Weighing Technology
GmbH
Registered Office:
Göttingen
Local Court of Registration:
Amtsgericht Göttingen
HRB 201458
Managing Directors:
Mario Hespe,
Carsten Wolter

Sartorius will continue to offer service and maintenance for Picus® and Picus® Nxt electronic pipettes for 5 years following discontinuation. After this 5-year period, normal calibration, upkeep service and spare parts will remain available, although availability of spare parts for repairs might be limited.

We continuously strive to implement solutions to our products and services to enable you to get your treatments to patients faster. We thank you for your trust and continued use of Sartorius products, and we look forward to assisting you in managing this change.

Sincerely,

DocuSigned by:
Fiona Coats
 Signer Name: Fiona Coats
 Signing Reason: I approve this document
 Signing Time: 2023-10-20 | 15:13:46 CEST
 9A53DBE65FC94520B4258D4D1B5CF7A7

DocuSigned by:
Susanne Gerighausen
 Name des Unterzeichners: Susanne Gerighausen
 Signiergrund: Ich genehmige dieses Dokument
 Signierzeit: 2023-10-20 | 15:09:18 MESZ
 165B30A4A29E4CD99853CE51CFCFDDBF

Fiona Coats
 Head of Marketing LPS
 Sartorius Corporate

Susanne Gerighausen
 Head of Quality Assurance & Control LPS
 Sartorius Lab Instruments GmbH & Co.KG

Note: List of affected manufacturing site(s) / manufacturing site(s) of impacted products
 -Sartorius Biohit Liquid Handling Oy, Helsinki, Finland
 -Sartorius Biohit Liquid Handling Oy, Kajaani, Finland

Appendix 1: List of affected products and alternative products

To be discontinued order number	Description	Alternative product order number	Description
	Picus® electronic pipettes		
735021	Picus® Electronic Pipette, 1-ch, 0.2 - 10 µL	LH-747021	Picus® 2 Electronic Pipette, 1-ch, 0.2 - 10 µL
735041	Picus® Electronic Pipette, 1-ch, 5 - 120 µL	LH-747041	Picus® 2 Electronic Pipette, 1-ch, 5 - 120 µL
735061	Picus® Electronic Pipette, 1-ch, 10 - 300 µL	LH-747061	Picus® 2 Electronic Pipette, 1-ch, 10 - 300 µL
735081	Picus® Electronic Pipette, 1-ch, 50 - 1,000 µL	LH-747081	Picus® 2 Electronic Pipette, 1-ch, 50 - 1,000 µL
735101	Picus® Electronic Pipette, 1-ch, 100 - 5,000 µL	LH-747101	Picus® 2 Electronic Pipette, 1-ch, 100 - 5,000 µL
735111	Picus® Electronic Pipette, 1-ch, 500 - 10,000 µL	LH-747111	Picus® 2 Electronic Pipette, 1-ch, 500 - 10,000 µL
735321	Picus® Electronic Pipette, 8-ch, 0.2 - 10 µL	LH-747321	Picus® 2 Electronic Pipette, 8-ch, 0.2 - 10 µL
735341	Picus® Electronic Pipette, 8-ch, 5 - 120 µL	LH-747341	Picus® 2 Electronic Pipette, 8-ch, 5 - 120 µL
735361	Picus® Electronic Pipette, 8-ch, 10 - 300 µL	LH-747361	Picus® 2 Electronic Pipette, 8-ch, 10 - 300 µL
735391	Picus® Electronic Pipette, 8-ch, 50 - 1,200 µL	LH-747391	Picus® 2 Electronic Pipette, 8-ch, 50 - 1,200 µL
735421	Picus® Electronic Pipette, 12-ch, 0.2 - 10 µL	LH-747421	Picus® 2 Electronic Pipette, 12-ch, 0.2 - 10 µL
735441	Picus® Electronic Pipette, 12-ch, 5 - 120 µL	LH-747441	Picus® 2 Electronic Pipette, 12-ch, 5 - 120 µL
735461	Picus® Electronic Pipette, 12-ch, 10 - 300 µL	LH-747461	Picus® 2 Electronic Pipette, 12-ch, 10 - 300 µL
735491	Picus® Electronic Pipette, 12-ch, 50 - 1,200 µL	LH-747491	Picus® 2 Electronic Pipette, 12-ch, 50 - 1,200 µL
	Picus® Nxt electronic pipettes		

To be discontinued order number	Description	Alternative product order number	Description
LH-745021	Picus® Nxt Electronic Pipette, 1-ch, 0.2 - 10 µL	LH-747021	Picus® 2 Electronic Pipette, 1-ch, 0.2 - 10 µL
LH-745041	Picus® Nxt Electronic Pipette, 1-ch, 5 - 120 µL	LH-747041	Picus® 2 Electronic Pipette, 1-ch, 5 - 120 µL
LH-745061	Picus® Nxt Electronic Pipette, 1-ch, 10 - 300 µL	LH-747061	Picus® 2 Electronic Pipette, 1-ch, 10 - 300 µL
LH-745081	Picus® Nxt Electronic Pipette, 1-ch, 50 - 1,000 µL	LH-747081	Picus® 2 Electronic Pipette, 1-ch, 50 - 1,000 µL
LH-745101	Picus® Nxt Electronic Pipette, 1-ch, 100 - 5,000 µL	LH-747101	Picus® 2 Electronic Pipette, 1-ch, 100 - 5,000 µL
LH-745111	Picus® Nxt Electronic Pipette, 1-ch, 500 - 10,000 µL	LH-747111	Picus® 2 Electronic Pipette, 1-ch, 500 - 10,000 µL
LH-745321	Picus® Nxt Electronic Pipette, 8-ch, 0.2 - 10 µL	LH-747321	Picus® 2 Electronic Pipette, 8-ch, 0.2 - 10 µL
LH-745341	Picus® Nxt Electronic Pipette, 8-ch, 5 - 120 µL	LH-747341	Picus® 2 Electronic Pipette, 8-ch, 5 - 120 µL
LH-745361	Picus® Nxt Electronic Pipette, 8-ch, 10 - 300 µL	LH-747361	Picus® 2 Electronic Pipette, 8-ch, 10 - 300 µL
LH-745391	Picus® Nxt Electronic Pipette, 8-ch, 50 - 1,200 µL	LH-747391	Picus® 2 Electronic Pipette, 8-ch, 50 - 1,200 µL
LH-745421	Picus® Nxt Electronic Pipette, 12-ch, 0.2 - 10 µL	LH-747421	Picus® 2 Electronic Pipette, 12-ch, 0.2 - 10 µL
LH-745441	Picus® Nxt Electronic Pipette, 12-ch, 5 - 120 µL	LH-747441	Picus® 2 Electronic Pipette, 12-ch, 5 - 120 µL
LH-745461	Picus® Nxt Electronic Pipette, 12-ch, 10 - 300 µL	LH-747461	Picus® 2 Electronic Pipette, 12-ch, 10 - 300 µL
LH-745491	Picus® Nxt Electronic Pipette, 12-ch, 50 - 1,200 µL	LH-747491	Picus® 2 Electronic Pipette, 12-ch, 50 - 1,200 µL