



Revvity, Inc.  
68 Elm St.  
Hopkinton, MA 01748  
USA  
Web: [www.revvity.com](http://www.revvity.com)

Technical  
Data Sheet

# LabChip® QC Test Results



### Lot Information

Lot Number: **F624B**  
Description: LabChip® 620N  
Test Date: **June 03, 2024**

Revvity hereby certifies that these products are manufactured in a controlled environment in accordance with established manufacturing guidelines, quality requirements and product specifications.

The above part number and associated lot number have passed in process and finished goods testing in accordance with the following testing and quality protocols.

| Parameter                                           | Specification                                         | Result |
|-----------------------------------------------------|-------------------------------------------------------|--------|
| Pressure Test                                       | Chips are free of clogs.                              | Pass   |
| Chip Interface                                      | Chips are recognized on LabChip® GX Touch Instrument. | Pass   |
| Autofocus                                           | Chips can autofocus on LabChip® GX Touch Instrument.  | Pass   |
| Upper Marker Migration Time                         | 49 to 62 seconds                                      | Pass   |
| Lower Marker-Upper Marker Migration Time Difference | $\geq 27$ seconds                                     | Pass   |
| Baseline Variability                                | $\leq 10\%$                                           | Pass   |
| Sensitivity                                         | $\geq 149$ RFU                                        | Pass   |
| Concentration Accuracy                              | $\pm 30\%$                                            | Pass   |
| Sizing Accuracy                                     | $\pm 10\%$                                            | Pass   |

It is hereby certified that the data and information in this document has been reviewed and is in conformance with applicable requirements and specifications for the part number and lot number above.