



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

Technical  
Data Sheet

# LabChip<sup>®</sup> QC Test Results



### Lot Information

Lot Number: **F090P**  
Description: LabChip<sup>®</sup> 563N  
Test Date: **February 16, 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 <sup>®</sup> GX Touch Instrument. | Pass   |
| Autofocus                   | Chips can autofocus on LabChip <sup>®</sup> GX Touch Instrument.  | Pass   |
| Lower Marker Migration Time | 11 to 13.5 seconds  | Pass   |
| Last Peak Migration Time    | 26 to 33 seconds  | Pass   |
| Last peak FWHM              | $\leq 0.22$   | Pass   |
| Concentration CV            | $\leq 25\%$   | Pass   |
| Peak Detect                 | All system and Ladder peaks detected                              | Pass   |
| Baseline                    | $\leq 250$ rfu  | 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.