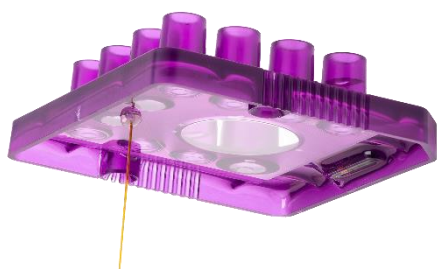




Revvity, Inc.  
68 Elm St.  
Hopkinton, MA 01748  
USA  
Web: [www.revvity.com](http://www.revvity.com)

Technical  
Data Sheet

# LabChip<sup>®</sup> QC Test Results



### Lot Information

Lot Number: **D392P**  
Description: LabChip<sup>®</sup> 563N  
Test Date: **September 21, 2023**

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.