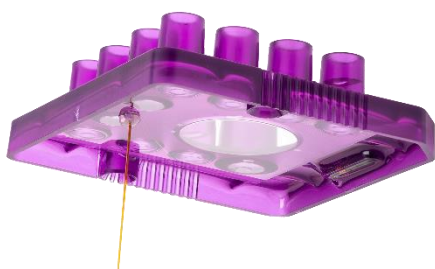


Technical
 Data Sheet

LabChip® QC Test Results



Lot Information

Lot Number: **F463B**
 Description: LabChip® 564N
 Test Date: **April 22, 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	20 to 25 seconds	Pass
Lower Marker-Upper Marker Migration Time Difference	>=8.5 seconds	Pass
Baseline Variability	<=9.5%	Pass
Concentration Accuracy	±30%	Pass
Sizing Accuracy	±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.