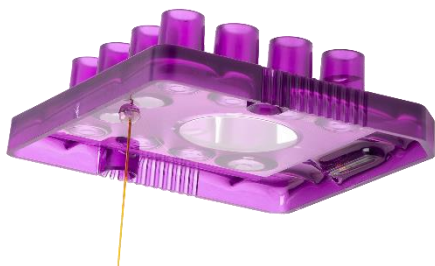




Revvity, Inc.
68 Elm St.
Hopkinton, MA 01748
USA
Web: www.revvity.com

Technical
Data Sheet

LabChip® QC Test Results



Lot Information

Lot Number: **F904P**
Description: LabChip® 584N
Test Date: **July 26, 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
Peak Detection	All system and ladder peaks are detected.	Pass
Last Peak Migration Time Average	22.5 to 27.5 seconds	Pass
Last peak Migration Time CV	$\leq 10\%$	Pass
Lower Marker Migration Time	17 to 20 seconds	Pass
FWHM	≤ 0.23	Pass
Area CV	$\leq 20\%$	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.