

Pipetting

Device qualification of the Finnpipette Novus electronic pipette (IQ/OQ/PQ)

Keywords

- Thermo Scientific Finnpipette Novus electronic pipette
- Device qualification
- Installation qualification (IQ)
- Operational qualification (OQ)
- Performance qualification (PQ)

Introduction

During device qualification, one verifies and documents whether the equipment, in this case a pipette, is reliable and capable of fulfilling the requirements of its intended use. All qualification activities are fully documented according to pertinent quality and regulatory requirements and signed by authorized personnel.

Installation qualification (IQ)

The main purpose of the IQ is to verify that the correct pipette has been delivered and that the electrical connections conform to the manufacturer's specifications and requirements.

IQ includes but is not limited to:

Verifications

- The catalog number on the label of the pipette shipping box corresponds to the purchase order.
- The pipette corresponds to the order specifications (volume range, number of channels, model).
- The catalog number stated on the Calibration Report corresponds to the purchase order.
- The serial number stated on the Calibration Report corresponds to the serial number on the pipette.
- The Instructions for Use corresponds to the purchased pipette model (either single channel or multichannel model as appropriate). The catalog number of the purchased model is stated in the Instructions for Use.
- The complete Thermo Scientific™ Finnpipette™



Novus™ product box contains the items stated in the Instructions for Use.

- The following environmental recommendations apply:
 - Operating temperature 15–35°C
 - Relative humidity during operation 20-85% and during storage 60 +/- 25%

Actions

- Confirm that the pipette charger is compatible with the local power supply and that the charger specifications correspond to the local power supply specifications.
- Fully charge the battery of the pipette, preferably for 2 hours, for the first charge.
- Confirm the existence of necessary SOPs (preventive maintenance, work instructions, etc.)
- Confirm that the required operator training has been conducted.

Operational qualification (OQ)

The OQ protocol verifies that the pipette is functioning in an acceptable manner. This includes testing the pipette to ensure that it is capable of operating in accordance with the specifications and requirements defined by the manufacturer.

OQ shall be performed after IQ and shall be repeated throughout the life of the pipette and after service. OQ must be included in the SOPs.

OQ includes but is not limited to:

- Ensuring the pipette is operating properly by checking basic operations and functions defined in the Instructions for Use, e.g. the function of the display, tip ejector and menu.
- Verifying pipette operation either by the end user or by an authorized party, e.g. the manufacturer. The Calibration Report provided with the pipette verifies the OQ performed by the manufacturer. User verification should be performed in a controlled environment using water, e.g. in accordance with ISO 8655. We recommend that end users perform the verifications themselves because pipetting results obtained by the user may differ from the values stated in the Calibration Report.
- Ensuring repeatable results, the pipette must be regularly serviced according to the Instructions for Use.

thermo scientific		PIPETTE CALIBRATION REPORT	
Product code:	4320000		
Description:	Finnpipette Nova 100-1000µl	Model ID:	
Pipette ID:	527873		
Tip:	FinnTip 1000		
No of channels:	1		
Manufacturer: Thermo Fisher Scientific, Järvenpää, Finland			
Tested by: 4327			
Date:	30.11.2021	Time:	11:45:55
Balance:	Ohaus N11105	Balance ID:	9120000
Environment data:	Ambio 2100	Env. data ID:	911000005
Air temperature:	22.85 °C	Water temperature:	20.82 °C
Humidity:	Moisture trap in use	Air pressure:	989.97Pa
Z-factor:	1.0032		
Volume:	100 µl	1000 µl	
1	100.0	1	1000.0
2	99.8	2	1000.0
3	100.1	3	999.9
4	99.9	4	999.9
5	100.0	5	999.9
Mean:	100.0 µl	999.9 µl	
s:	0.4 µl	0.4 µl	
k:	0.7 µl	0.9 µl	
AccL:	0.24	0.04	
CV%	0.12	0.09	
	PASS500	PASS500	
Manufacturer specifications:			
Min volume:	10 µl	Max volume:	1000 µl
A:	0.4 µl		0 µl
k:	0.6 µl		2 µl
Acc%	0.20		0.05
CV%	0.80		0.20
Manufacturer specifications meet the requirements of ISO 8655			

Example of Calibration Report.

Performance qualification (PQ)

The purpose of the PQ is to verify that the pipette is capable of properly accomplishing the task for its intended use. PQ relates to the daily use of the pipette and is designed to measure routine performance.

Note: Prior to shipment, the pipettes are tested under factory conditions. However, performance in a laboratory may vary due to temperature, pressure, humidity, operator, liquid, type of tip, etc. This should be taken into account when verification and validation tests are performed.

PQ includes but is not limited to:

- Determination by the user of the acceptance criteria required by the most critical pipette and tip application. Typically, this means setting acceptance limits for pipetting accuracy and precision and verifying that the pipette and tip fulfill these criteria.

Because the pipette typically is part of a broad application comprised of several instruments, system validation is recommended. This should be done after PQ has been performed for each instrument. In practice, system validation often requires the use of proven calibrators, samples or reference liquids.

