



### CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** MESO® QuickPlex Q 60MM Performance Qualification Kit

**Catalog Number:** RA2QQ-1

**Lot Number:** PQ500010

**Expiration Date:** 30 NOV 2022

**Storage Temperature:** Room Temperature

Components	Catalog Number	Lot Number	Expiration Date
MULTI-ARRAY 96-Well QUICKPLEX Plate	L55XA	2MK0HA	30 NOV 2022
PQ Control 1	N/A	Y0260166	31 MAY 2023
PQ Negative Control	R92TG	Y0600015	31 MAY 2023

Test	QC Obtained Mean Signal	Signal Specifications	
		Minimum Value	Maximum Value
PQ Control 1	14791	13312	16270
PQ Negative Control	72	37	107
Dark Noise	0	-16	16
Test	%CV and SD Specification		
PQ Control 1	0-6.0%		
PQ Negative Control	0-25		
Dark Noise	0-16		

This product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
<b>Review/Approval</b>	Karen Hamilla	Quality		03 MAR 2021