



### CERTIFICATE OF ANALYSIS

Page 1 of 1

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

**Catalog Number:** RA2QQ-1

**Lot Number:** PQ500008

**Expiration Date:** 30 SEP 2021

**Storage Temperature:** Room Temperature

Components	Catalog Number	Lot Number	Expiration Date
MULTI-ARRAY 96-Well QUICKPLEX Plate	L55XA	2MJ0FA	30 SEP 2021
PQ Control 1	N/A	Y0260165	31 DEC 2022
PQ Negative Control	R92TG	N/A	N/A

Test	QC Obtained Mean Signal	Signal Specifications	
		Minimum Value	Maximum Value
PQ Control 1	14308	12877	15739
PQ Negative Control	73	38	108
Dark Noise	1	-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		13 JAN 2020