



CERTIFICATE OF ANALYSIS

Page 1 of 1

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

Catalog Number: RA2QQ-1

Lot Number: PQ500016

Expiration Date: 30 NOV 2024

Storage Temperature: Room Temperature

Specifications:

Specification	Minimum	Maximum
PQ Control 1 Mean Signal	15398	18820
PQ Control 1 %CV	0	6.0%
PQ Negative Control Mean Signal	48	118
PQ Negative Control Standard Deviation	0	25
Dark Noise Mean Signal	-16	16
Dark Noise Standard Deviation	0	16

MSD Testing Results:

Metric	Observed
PQ Control 1 Median of Mean Signal	17109
PQ Negative Control Median of Mean Signal	83
Dark Noise Median of Mean Signal	-4

Lot-specific Components:

Description	Catalog Number	Lot Number	Expiration Date
QuickPlex® 96-Well Plate	L55XA	2MM0PA	30 NOV 2024
PQ Control 1	N/A	Y0260176	31 AUG 2025
PQ Negative Control	N/A	Y0600037	30 JUN 2025

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		29 JUN 2023