



CERTIFICATE OF ANALYSIS

Product Description: MESO® QuickPlex Q 60MM Performance Qualification Kit
Catalog Number: RA2QQ-1
Lot Number: PQ500017
Expiration Date: 31 AUG 2025
Storage Temperature: Room Temperature

Lot-Specific Components:

Description	Lot Number	Expiration Date
QuickPlex® 96-Well Plate	2MN0GA	31 AUG 2025
PQ Control 1	Y0260177	31 OCT 2025
PQ Negative Control	Y0600044	31 AUG 2026

MSD Testing Results:

Metric	Observed
PQ Control 1 Median of Mean Signal	14581
PQ Negative Control Median of Mean Signal	75
Dark Noise Median of Mean Signal	-1

Specifications:

Specification	Minimum	Maximum
PQ Control 1 Mean Signal	13123	16039
PQ Control 1 %CV	0	6%
PQ Negative Control Mean Signal	40	110
PQ Negative Control Standard Deviation	0	25
Dark Noise Mean Signal	-16	16
Dark Noise Standard Deviation	0	16

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		28 FEB 2024