

**CERTIFICATE OF ANALYSIS**

Product Description: MESO QuickPlex Q 60MM Performance Qualification Kit
Catalog Number: RA2QQ-1
Lot Number: PQ500023
Expiration Date: 30 APR 2027
Storage Temperature: Room Temperature

Lot-Specific Components:

Description	Lot Number	Expiration Date
QuickPlex® 96-Well Plate	2MP0DA	30 APR 2027
PQ Control 1	Y0260183	31 DEC 2027
PQ Negative Control	Y0600052	30 JUN 2027

MSD Testing Results:

Metric	Observed
PQ Control 1 Median of Mean Signal	15997
PQ Negative Control Median of Mean Signal	84
Dark Noise Median of Mean Signal	-1

Specifications:

Specification	Minimum	Maximum
PQ Control 1 Mean Signal	14398	17597
PQ Control 1 %CV	0	6%
PQ Negative Control Mean Signal	49	119
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		30 JUN 2026