



CERTIFICATE OF ANALYSIS

Page 1 of 1

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

Catalog Number: RA2QQ-1

Lot Number: PQ500011

Expiration Date: 30 APR 2023

Storage Temperature: Room Temperature

Specifications:

Specification	Minimum	Maximum
PQ Control 1 Mean Signal	13983	17090
PQ Control 1 %CV	0	6.0%
PQ Negative Control Mean Signal	47	117
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	15537
PQ Negative Control Median of Mean Signal	82
Dark Noise Median of Mean Signal	1

Lot-specific Components:

Description	Catalog Number	Lot Number	Expiration Date
QuickPlex® 96-Well Plate	L55XA	2ML06A	30 APR 2023
PQ Control 1	N/A	Y0260166	31 MAY 2023
PQ Negative Control	N/A	Y0600024	30 JUN 2024

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

	Name	Function	Signature	Date
Review/Approval	Carissa B. Hubbard	Quality		20 AUG 2021