



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: S 600/600MM and SQ 120/120MM Performance Qualification Kit

Catalog Number: R31QQ-3

Lot Number: PQ300154

Expiration Date: 30 SEP 2024

Storage Temperature: Room Temperature

Specifications:

Specification	Minimum	Maximum
Free Tag Mean Signal	12000	18000
Free Tag %CV	0	6.0%
PQ Low Control Mean Signal	25.0	100
PQ Low Control Standard Deviation	0	25
Dark Noise Mean Signal	-16	16
Dark Noise Standard Deviation	0	16

MSD Testing Results:

Metric	Observed
Free Tag Mean Signal	15131
PQ Low Control Mean Signal	64
Dark Noise Mean Signal	2

Lot-specific Components:

Description	Catalog Number	Lot Number	Expiration Date
MULTI-ARRAY® 96-well Plate	L15XA	21M4FA	30 SEP 2024
MSD Free Tag ECL 15 000	N/A	Y0260177	31 OCT 2025
PQ Low Control	R92TE	Y0110095	30 NOV 2025

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		23 JAN 2023