



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: PQ300144

Expiration Date: 31 AUG 2023

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	15995
PQ Low Control Mean Signal	68
Dark Noise Mean Signal	-2

Lot-specific Components:

Description	Catalog Number	Lot Number	Expiration Date
MULTI-ARRAY® 96-well Plate	L15XA	21L3UA	31 AUG 2023
MSD Free Tag ECL 15 000	N/A	Y0260171	31 MAY 2024
PQ Low Control	R92TE	Y0110089	30 JUN 2024

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		29 SEP 2021