



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: PQ300140

Expiration Date: 31 JUL 2022

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	15778
PQ Low Control Mean Signal	70.0
Dark Noise Mean Signal	3.0

Lot-specific Components:

Description	Catalog Number	Lot Number	Expiration Date
MULTI-ARRAY® 96-well Plate	L15XA	21K2WA	31 JUL 2022
MSD Free Tag ECL 15 000	N/A	Y0260167	31 AUG 2023
PQ Low Control	R92TE	Y0110085	30 NOV 2023

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		23 DEC 2020