



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: PQ300141

Expiration Date: 28 FEB 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	15486
PQ Low Control Mean Signal	75.0
Dark Noise Mean Signal	1.0

Lot-specific Components:

Description	Catalog Number	Lot Number	Expiration Date
MULTI-ARRAY® 96-well Plate	L15XA	21L0CA	28 FEB 2023
MSD Free Tag ECL 15 000	N/A	Y0260169	31 DEC 2023
PQ Low Control	R92TE	Y0110086	31 JAN 2024

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		31 MAR 2021