



## 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:** PQ300149

**Expiration Date:** 31 MAR 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	15346
PQ Low Control Mean Signal	66
Dark Noise Mean Signal	0

### Lot-specific Components:

Description	Catalog Number	Lot Number	Expiration Date
MULTI-ARRAY® 96-well Plate	L15XA	21M13A	31 MAR 2024
MSD Free Tag ECL 15 000	N/A	Y0260173	31 DEC 2024
PQ Low Control	R92TE	Y0110092	31 JAN 2025

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		02 JUN 2022