



## 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:** PQ300155

**Expiration Date:** 31 JAN 2025

**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	14214
PQ Low Control Mean Signal	59
Dark Noise Mean Signal	2

### Lot-specific Components:

Description	Catalog Number	Lot Number	31 JAN 2025
MULTI-ARRAY® 96-well Plate	L15XA	21N0CA	31 JAN 2025
MSD Free Tag ECL 15 000	N/A	Y0260177	31 OCT 2025
PQ Low Control	R92TE	Y0110094	31 MAY 2025

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		28 FEB 2023