



CERTIFICATE OF ANALYSIS

**Product Description:** S 600/600MM and SQ 120/120MM Performance Qualification Kit  
**Catalog Number:** R31QQ-3  
**Lot Number:** PQ300167  
**Expiration Date:** 31 MAY 2026  
**Storage Temperature:** Room Temperature

**Lot-Specific Components:**

Description	Lot Number	Expiration Date
MULTI-ARRAY® 96-Well Plate	21O35A	31 MAY 2026
MSD Free Tag ECL 15000	Y0260181	31 MAY 2027
PQ Low Control	Y0320016	31 MAY 2027

**MSD Testing Results:**

Metric	Observed
Free Tag Mean Signal	15175
PQ Low Control Mean Signal	73
Dark Noise Mean Signal	1

**Specifications:**

Specification	Minimum	Maximum
Free Tag Mean Signal	12000	18000
Free Tag %CV	0	6%
PQ Low Control Mean Signal	25	100
PQ Low Control Standard Deviation	0	25
Dark Noise Mean Signal	-16	16
Dark Noise Standard Deviation	0	16

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		21 AUG 2024