



### CERTIFICATE OF ANALYSIS

**Product Description:** S 600/600MM and SQ 120/120MM Performance Qualification Kit  
**Catalog Number:** R31QQ-3  
**Lot Number:** PQ300182  
**Expiration Date:** 31 OCT 2027  
**Storage Temperature:** Room Temperature

**Lot-Specific Components:**

Description	Lot Number	Expiration Date
96-Well 1-Spot SECTOR Plate	21P5OA	31 OCT 2027
MSD Free Tag ECL 15000	Y0260186	31 JAN 2029
PQ Low Control	Y0320018	30 SEP 2028

**MSD Testing Results:**

Metric	Observed
Free Tag Mean Signal	14360
PQ Low Control Mean Signal	56
Dark Noise Mean Signal	4

**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
<b>Review/Approval</b>	Karen Hamilla	Quality		17 MAR 2026