

**CERTIFICATE OF ANALYSIS**

Product Description: S 600/600MM and SQ 120/120MM Performance Qualification Kit
Catalog Number: R31QQ-3
Lot Number: PQ300171
Expiration Date: 30 NOV 2026
Storage Temperature: Room Temperature

Lot-Specific Components:

Description	Lot Number	Expiration Date
MULTI-ARRAY® 96-Well Plate	21O77A	30 NOV 2026
MSD Free Tag ECL 15000	Y0260182	31 AUG 2027
PQ Low Control	Y0320017	31 DEC 2027

MSD Testing Results:

Metric	Observed
Free Tag Mean Signal	14449
PQ Low Control Mean Signal	65
Dark Noise Mean Signal	3

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		14 MAR 2025