



CERTIFICATE OF ANALYSIS

Product Description: SECTOR PR2 Model 1800 Performance Qualification Kit
Catalog Number: R61QQ-2
Lot Number: PQ400058
Expiration Date: 30 APR 2026
Storage Temperature: Room Temperature

Lot-Specific Components:

Description	Lot Number	Expiration Date
Uncoated Test Plate, 9 Spot 96-Well PR2 Plate	1VO0DA	30 APR 2026
MSD Free Tag ECL 15000	Y0260181	31 MAY 2027

MSD Testing Results:

Metric	Observed
Free Tag Mean Signal	11386

Specifications:

Specification	Minimum	Maximum
Free Tag Mean Signal	10000	20000
Free Tag %CV	0	15%

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		24 JUL 2024