



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: SECTOR® PR2 Model 1800 Performance Qualification Kit

Catalog Number: R61QQ-2

Lot Number: PQ400053

Expiration Date: 30 SEP 2023

Storage Temperature: Room Temperature

Specifications:

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

MSD Testing Results:

Metric	Observed
Free Tag Mean Signal	13982

Lot-specific Components:

Description	Catalog Number	Lot Number	Expiration Date
Uncoated Test Plate, 9 Spot 96-Well PR2 Plate	2072-0001	1VL0HA	30 SEP 2023
MSD Free Tag ECL 15000	N/A	Y0260172	31 AUG 2024

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		18 FEB 2022