



### CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** SECTOR® PR2 Model 1800 Performance Qualification Kit

**Catalog Number:** R61QQ-2

**Lot Number:** PQ400054

**Expiration Date:** 30 JUN 2024

**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	14330

#### Lot-specific Components:

Description	Catalog Number	Lot Number	Expiration Date
Uncoated Test Plate, 9 Spot 96-Well PR2 Plate	2072-0001	1VM06A	30 JUN 2024
MSD Free Tag ECL 15 000	N/A	Y0260174	31 MAR 2025

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		18 AUG 2022