



### CERTIFICATE OF ANALYSIS

Page 1 of 1

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

**Catalog Number:** R61QQ-2

**Lot Number:** PQ400046

**Expiration Date:** 31 AUG 2021

**Storage Temperature:** Room Temperature

Components	Catalog Number	Lot Number	Expiration Date
Uncoated Test Plate, 9 Spot 96-Well PR2 Plate	2072-0001	1VJ1JA	31 AUG 2021
MSD Free Tag ECL 15 000	N/A	Y0260164	30 SEP 2022

Test	QC Obtained Mean Signal	Signal Specifications	
		Minimal Value	Maximal Value
MSD Free Tag ECL 15000	13573	10 000	20 000
	%CV Specification		
	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 OCT 2019