



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: SECTOR® PR2 Model 1800 Performance Qualification Kit

Catalog Number: R61QQ-2

Lot Number: PQ400048

Expiration Date: 31 JUL 2021

Storage Temperature: Room Temperature

Components	Catalog Number	Lot Number	Expiration Date
Uncoated Test Plate, 9 Spot 96-Well PR2 Plate	2072-0001	1VJ14A	31 JUL 2021
MSD Free Tag ECL 15 000	N/A	Y0260166	31 MAY 2023

Test	QC Obtained Mean Signal	Signal Specifications	
		Minimal Value	Maximal Value
MSD Free Tag ECL 15000	16133	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		21 SEP 2020