



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: SECTOR[®] PR2 Model 1800 Performance Qualification Kit 2

Catalog Number: R81QQ-1

Lot Number: PQ600001

Expiration Date: 30 NOV 2022

Storage Temperature: Room Temperature

Kit Components:

Description	Lot Number	Expiration Date
PR2 PQ Test Plate	2MK00A	30 NOV 2022
PR2 PQ Control 1	Y0810001	30 JUN 2023

Functional Testing Results:

Metric	Value
Mean Signal	14727

PQ Kit Specifications:

Specification	Value
Maximum Signal	17672
Minimum Signal	11782
Maximum % CV	10%

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		16 MAR 2021