



CERTIFICATE OF ANALYSIS

Page 1 of 1

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

Catalog Number: R81QQ-1

Lot Number: PQ600002

Expiration Date: 31 DEC 2024

Storage Temperature: Room Temperature

Kit Components:

Description	Lot Number	Expiration Date
PR2 PQ Test Plate	2MM0TA	31 DEC 2024
PR2 PQ Control 1	Y0810002	31 JUL 2025

Functional Testing Results:

Metric	Value
Mean Signal	10693

PQ Kit Specifications:

Specification	Value
Maximum Signal	12832
Minimum Signal	8554
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		07 MAR 2023