



CERTIFICATE OF ANALYSIS

Product Description: SECTOR® PR2 Model 1800 Performance Qualification Kit 2
Catalog Number: R81QQ-1
Lot Number: PQ600005
Expiration Date: 31 AUG 2025
Storage Temperature: Room Temperature

Kit Components:

Description	Lot Number	Expiration Date
PR2 PQ Test Plate	2MN0IA	31 AUG 2025
PR2 PQ Control 1	Y0810006	31 OCT 2026

Functional Testing Results:

Metric	Value
Mean Signal	14180

PQ Kit Specifications:

Specification	Value
Maximum Signal	17016
Minimum Signal	11344
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		17 JUN 2024