



CERTIFICATE OF ANALYSIS

COA – 04409

Product Description: MSD® Performance Qualification Kit, PR

Catalog #/Size: R21QQ-2

Lot Number: PQ100019

Expiration Date: 30 APR 2015

Statement: This product was manufactured and tested according to approved MSD documents. This lot meets MSD’s specifications.

SECTOR PR2 Plate Lot Number: 01D02A
SECTOR PR2 Plate Catalog Number: L13XA
SECTOR PR2 Plate Expiration Date: 30 APR 2015
Free Tag 15 000 Lot Number: Y0260143
Free Tag 15 000 Part Number: N/A
Free Tag 15 000 Expiration Date: 31 AUG 2016
Read Buffer T(1x) Lot Number: Y0110049
Read Buffer T(1x) Catalog Number: R92TA
Read Buffer T(1x) Expiration Date: 31 MAY 2016

Specifications:

Table with 3 columns: Parameter, Max, Min. Rows include Free Tag Mean, Free Tag CV %, Read Buffer T(1x) Mean, Read Buffer T(1x) SD, Dark Signal Mean, and Dark Noise (SD).

MSD QC Results: Using instrument: PR400 S/N: 1150040915035
Free TAG Mean: 15 193
Read Buffer T(1X) Mean: 66
Dark Signal Mean: 0

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

Table with 5 columns: Role, Name, Function, Signature, Date. Rows for Preparer (Rolanda Spratley), Review/Approval (Micheal Eager), and Review/Approval (Richard Mann).