



CERTIFICATE OF ANALYSIS

COA – 04273

Product Description: MSD® Performance Qualification Kit, Sector Imager

Catalog #/Size: R31QQ-3

Lot Number: PQ300039

Expiration Date: 31 MAR 2016

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

- SECTOR Imager Plate Lot Number: 21E0HA
SECTOR Imager Plate Catalog Number: L15XA
SECTOR Imager Plate Expiration Date: 31 MAR 2016
Free Tag 15,000 Lot Number: Y0260145
Free Tag 15,000 Catalog Number: N/A
Free Tag 15,000 Expiration Date: 31 JAN 2017
PQ Low Control Lot Number: Y0320002
PQ Low Control Catalog Number: R92TE
PQ Low Control Expiration Date: 31 JAN 2017

Specifications:

Table with 3 columns: Parameter, Max, Min. Rows include Free Tag Mean, Free Tag CV, PQ Low Control Mean, PQ Low Control Mean SD, Dark Signal Mean, and Dark Noise (SD).

MSD QC Results: Using instrument SI: 6000 S/N: 1200050211221
Free TAG Mean: 16 535
PQ Low Control Mean: 80
Dark Signal Mean: -1

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 include Preparer (Rolanda Spratley), Review/Approval (Michael Eager), and Review/Approval (Richard Mann).