



CERTIFICATE OF ANALYSIS

COA - 07209

Product Description: MSD® Bioprocess Duplex Kit

Catalog #/Size: K15149F-1 / 1 Plate; K15149F-2 / 5 Plates; K15149F-3 / 20 Plates

Lot Number: K0050714

Expiration Date: N/A (See expiration date on specific components below)

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

Kit Contents:

- Plate Lot Number: Z0055397
Plate Lot Catalog Number: N45149B
Plate Lot Expiration Date: 31 JAN 2018
Antibody blend Lot Number: D0050641
Antibody blend Catalog Number: D20HM
Antibody blend Expiration Date: 31 JAN 2018
Antibody blend Lot Number: D0050640
Antibody blend Catalog Number: D20HN
Antibody blend Expiration Date: 31 JAN 2018
Calibrator blend Lot Number: A0020179
Calibrator blend Catalog Number: C00HM
Calibrator blend Expiration Date: 31 JAN 2018
Calibrator blend Lot Number: A0020178
Calibrator blend Catalog Number: C00HN
Calibrator blend Expiration Date: 31 JAN 2018

Specifications:

- Release method: Coating uniformity
Mean Intraplate CV: ≤ 15%
Mean Interplate CV: ≤ 18%
No more than one plate with CV > 18%
No plate with CV > 20%

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

Table with 5 columns: Name, Function, Signature, Date. Rows include Preparer (Xiomara Talavera) and Review/Approval (Vanessa Castleberry).