



CERTIFICATE OF ANALYSIS

COA - 07269

Product Description: MSD® Human IFN-alpha2a Ultra-Sensitive Kit
Catalog #/Size: K151ACC-1 / 1 Plate; K151ACC -2 / 5 Plates; K151ACC -4 / 25 Plates
Lot Number: K0035207
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: Z0046150
Plate Lot Catalog Number: N45032A
Plate Lot Expiration Date: 30 NOV 2018
Antibody blend Lot Number: D0033664
Antibody blend Catalog Number: D21AC
Antibody blend Expiration Date: 30 NOV 2018
Antibody source (host/species): Mouse
Calibrator blend Lot Number: A0032247
Calibrator blend Catalog Number: C0032
Calibrator blend Expiration Date: 30 NOV 2018
Calibrator source (host/species): E. coli

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 (Dawn Stollar).