



CERTIFICATE OF ANALYSIS

COA - 07051

Product Description: MSD® Human IFN-alpha2a Tissue Culture Kit
Catalog #/Size: K151ACB -1 / 1 Plate; K151ACB-2 / 5 Plates; K151ACB-4 / 25 Plates
Lot Number: K0035134
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: Z0046063
Plate Lot Catalog Number: N45032B
Plate Lot Expiration Date: 31 AUG 2018
Antibody blend Lot Number: D0033610
Antibody blend Catalog Number: D21AC
Antibody blend Expiration Date: 31 AUG 2018
Antibody source (host/species): Mouse
Calibrator blend Lot Number: A0032225
Calibrator blend Catalog Number: C0032
Calibrator blend Expiration Date: 31 AUG 2018
Calibrator source (host/species): h1L-12(p40) - Sf 21
hI-TAC; hIFNalpha -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 (Rolanda Spratley) and Review/Approval (Dawn Stollar).