



CERTIFICATE OF ANALYSIS

COA - 06219

Product Description: MSD[®] Human IFN-gamma Tissue Culture Kit
Catalog #/Size: K151AEB-1 / 1 Plate; K151AEB-2 / 5 Plates; K151AEB-4 / 25 Plates
Lot Number: K0035000
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: Z0045820
 Plate Lot Catalog Number: L451AEB
 Plate Lot Expiration Date: 31 JAN 2018
 Antibody blend Lot Number: D0033491
 Antibody blend Catalog Number: D21AE
 Antibody blend Expiration Date: 31 JAN 2018
 Antibody source (host/species): Mouse
 Calibrator blend Lot Number: A0032157
 Calibrator blend Catalog Number: C01XX
 Calibrator blend Expiration Date: 28 FEB 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.

	Name	Function	Signature	Date
Preparer	Rolanda Spratley	Biologics/Chemistry		18 MAR 2016
Review/Approval	Dawn Stollar	Quality Assurance		22 MAR 2016