



CERTIFICATE OF ANALYSIS

COA - 07193

Product Description: MSD[®] Human IL-10 Tissue Culture Kit

Catalog #/Size: K151AOB-1 / 1 Plate; K151AOB-2 / 5 Plates; K151AOB-4 / 25 Plates

Lot Number: K0035174

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: Z0046123

Plate Lot Catalog Number: N45015B

Plate Lot Expiration Date: 30 SEP 2018

Antibody blend Lot Number: D0033641

Antibody blend Catalog Number: D21AO

Antibody blend Expiration Date: 30 SEP 2018

Antibody source (host/species): Rat

Calibrator blend Lot Number: A0032242

Calibrator blend Catalog Number: C0015

Calibrator blend Expiration Date: 30 SEP 2018

Calibrator source (host/species): Sf 21

Specifications:

Release method: Coating uniformity

Mean Intraplate CV: $\leq 15\%$

Mean Interplate CV: $\leq 18\%$

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	Xiomara Talavera	Quality Control		05 AUG 2016
Review/Approval	Vanessa Castleberry	Quality Assurance		08 AUG 2016