



CERTIFICATE OF ANALYSIS

COA - 06377

Product Description: MSD[®] Human Eotaxin-2 Kit

Catalog #/Size: K151RKD-1 / 1 Plate; K151RKD-2 / 5 Plates; K151RKD-4 / 25 Plates

Lot Number: K00E0281

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: Z00E0229

Plate Lot Catalog Number: N451RKA

Plate Lot Expiration Date: 30 NOV 2018

Antibody blend Lot Number: D00E0262

Antibody blend Catalog Number: D21RK

Antibody blend Expiration Date: 30 NOV 2018

Calibrator blend Lot Number: A00E0141

Calibrator blend Catalog Number: C01RK

Calibrator blend Expiration Date: 31 DEC 2018

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	<i>Xiomara Talavera</i>	08 SEP 2016
Review/Approval	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	08 SEP 2016