



CERTIFICATE OF ANALYSIS

COA - 07902

Product Description: MSD[®] Human IL-18 Kit

Catalog #/Size: K151MCD - Series

Lot Number: K00E0282

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: Z00E0228

Plate Lot Catalog Number: N451MCA

Plate Lot Expiration Date: 30 NOV 2018

Antibody blend Lot Number: D00E0263

Antibody blend Catalog Number: D21MC

Antibody blend Expiration Date: 30 NOV 2018

Calibrator blend Lot Number: A00E0142

Calibrator blend Catalog Number: C01MC

Calibrator blend Expiration Date: 30 NOV 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		18 FEB 2017
Review/Approval	Vanessa Castleberry	Quality Assurance		21 FEB 2017