



## CERTIFICATE OF ANALYSIS

COA - 06998

**Product Description:** MSD<sup>®</sup> Human IL-18 Kit

**Catalog #/Size:** K151MCD-1 / 1 Plate; K151MCD-2 / 5 Plates; K151MCD-4 / 25 Plates

**Lot Number:** K00E0204

**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: Z00E0177

Plate Lot Catalog Number: N451MCA

Plate Lot Expiration Date: 31 OCT 2017

Antibody blend Lot Number: D00E0197

Antibody blend Catalog Number: D21MC

Antibody blend Expiration Date: 30 SEP 2017

Calibrator blend Lot Number: A00E0105

Calibrator blend Catalog Number: C01MC

Calibrator blend Expiration Date: 30 SEP 2017

### 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
<b>Preparer</b>	Rolanda Spratley	Biologics/Chemistry		08 JUN 2016
<b>Review/Approval</b>	Vanessa Castleberry	Quality Assurance		08 JUN 2016