



### CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** MSD® Human MMP 2-Plex Ultra-Sensitive Kit

**Catalog Number:** K15033C-1 / 1 Plate; K15033C-2 / 5 Plates; K15033C-4 / 25 Plates

**Lot Number:** K0035105

**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					
<b>Lot</b>	Z0046028	<b>Catalog #</b>	N45033A	<b>Expiration Date</b>	31 JUL 2018

Antibody Blend					
<b>Lot</b>	D0033586	<b>Catalog #</b>	D2033	<b>Expiration Date</b>	31 JUL 2018
<b>Spot #</b>	<b>Description</b>		<b>Host / Species</b>		
1	hMMP-2		Goat		
2	hMMP-10		Goat		

Calibrator Blend					
<b>Lot</b>	A0032211	<b>Catalog #</b>	C0033	<b>Expiration Date</b>	31 AUG 2018
<b>Spot #</b>	<b>Description</b>		<b>Host / Species</b>		
1	hMMP-10		NS0		
2	hMMP-2		NS0		

Specifications:

Release method:

Coating Uniformity	Specifications
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
<b>Preparer</b>	Xiomara Talavera	Quality Control		04 NOV 2016
<b>Review/Approval</b>	Dawn Stollar	Quality Assurance		04 NOV 2016