



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: MSD® Human MMP 3-Plex Ultra-Sensitive Kit

Catalog Number: K15034C-1 / 1 Plate; K15034C-2 / 5 Plates; K15034C-4 / 25 Plates

Lot Number: K0035162

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	Z0046090	Catalog #	N45034A	Expiration Date	31 DEC 2018

Antibody Blend					
Lot	D0033631	Catalog #	D2034	Expiration Date	30 NOV 2018
Spot #	Description		Host / Species		
1	hMMP-3		Goat		
2	hMMP-9		Goat		
3	hMMP-1		Goat		

Calibrator Blend					
Lot	A0032248	Catalog #	C0034	Expiration Date	31 DEC 2018
Spot #	Description		Host / Species		
1	hMMP-1		NS0		
2	hMMP-3		NS0		
3	hMMP-9		CHO		

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
Preparer	Xiomara Talavera	Quality Control	<i>Xiomara Talavera</i>	04 NOV 2016
Review/Approval	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	04 NOV 2016