



### CERTIFICATE OF ANALYSIS

Page 1 of 1

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

**Catalog Number:** K15034C-Series

**Lot Number:** K0035731

**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>	Z0046838	<b>Catalog #</b>	N45034A	<b>Expiration Date:</b>	30 NOV 2020

Antibody Blend	
<b>Lot Number:</b>	D0034069
<b>Catalog Number:</b>	D2034
<b>Expiration Date:</b>	31 MAR 2022
<b>Storage Temperature:</b>	2-8°C
<b>Spot #</b>	<b>Description</b>
1	Human MMP-1
2	Human MMP-3
3	Human MMP-9

Calibrator Blend	
<b>Lot Number:</b>	A0032397
<b>Catalog Number:</b>	C0034
<b>Expiration Date:</b>	30 NOV 2020
<b>Storage Temperature:</b>	≤ -70°C
<b>Spot #</b>	<b>Description</b>

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>Review/Approval</b>	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	21 DEC 2018