



CERTIFICATE OF ANALYSIS

Page 1 of 1

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

Catalog Number: K151FZC-Series

Lot Number: K0035571

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 #	Z0046628	Catalog #	N45034A	Expiration Date:	28 FEB 2020

Antibody Blend	
Lot Number:	D0033936
Catalog Number:	D2034
Expiration Date:	29 FEB 2020
Storage Temperature:	2-8°C
Spot #	Description
1	Human MMP-1
2	Human MMP-3
3	Human MMP-9

Calibrator Blend	
Lot Number:	A0032333
Catalog Number:	C0034
Expiration Date:	28 FEB 2020
Storage Temperature:	≤ -70°C
Spot #	Description
1	Human MMP 3-Plex

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
Review/Approval	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	21 NOV 2017