



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: MSD® Isotyping Panel 1 (Human/NHP) Kit

Catalog Number: K15203D - Series

Lot Number: K00E0414

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 #	Z00E0324	Catalog #	N45203A	Expiration Date:	31 JUL 2020

Antibody Blend	
LotNumber:	D00E0375
Catalog Number:	D20JJ
Expiration Date:	30 NOV 2021
Storage Temperature:	2-8°C

Calibrator Blend	
Lot Number:	A00E0202
Catalog Number:	C0203
Expiration Date:	31 AUG 2020
Storage Temperature:	≤ -70°C

Antibody Blend	
LotNumber:	D00E0387
Catalog Number:	D20JL
Expiration Date:	31 JUL 2021
Storage Temperature:	2-8°C

Antibody Blend	
LotNumber:	D00E0388
Catalog Number:	D20JP
Expiration Date:	31 JUL 2020
Storage Temperature:	2-8°C

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>	26 FEB 2019