



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: MSD® Human YKL-40 Kit

Catalog Number: K151NHD-Series

Lot Number: K00E0361

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 #	Z00E0283	Catalog #	N451NHA	Expiration Date:	31 DEC 2019

Antibody Blend	
LotNumber:	D00E0336
Catalog Number:	D21NH
Expiration Date:	31 DEC 2019
Storage Temperature:	2-8°C

Calibrator Blend	
Lot Number:	A00E0181
Catalog Number:	C01NH
Expiration Date:	31 DEC 2019
Storage Temperature:	≤ -70°C

Specifications:

Release method:

Coating Uniformity	Specifications
Mean Intraplate CV	≤ 15%
No more than one plate with CV	> 18%
No plate with CV	> 25%

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>	27 AUG 2018