

CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: MSD® Human YKL-40 Kit

Catalog Number: K151NHD-Series

Lot Number: K00E0155

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 #	Z00E0142	Catalog #	N451NHA	Expiration Date:	30 NOV 2016

Antibody Blend		
LotNumber:	D00E0146	
Catalog Number:	D21NH	
Expiration Date:	31 DEC 2016	
Storage Temperature:	2-8°C	

Calibrator Blend			
Lot Number:	A00E0084		
Catalog Number:	C01NH		
Expiration Date:	30 NOV 2016		
Storage Temperature:	≤ -70°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	Vanessa Castleberry	Quality Assurance	Vanessa Castlebury	18MAY2017

COA-8271-A QA-FM-153-A