



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: MSD® Human Angiopoietin-1 Kit

Catalog Number: K151LPD-Series

Lot Number: K00E0362

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 #	Z00E0282	Catalog #	N451LPA	Expiration Date:	31 DEC 2019

Antibody Blend	
LotNumber:	D00E0337
Catalog Number:	D21LP
Expiration Date:	31 DEC 2019
Storage Temperature:	2-8°C

Calibrator Blend	
Lot Number:	A00E0182
Catalog Number:	C01LP
Expiration Date:	31 DEC 2019
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	<i>Vanessa Castleberry</i>	13SEP2017