



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Human Angiopoietin-2 Kit
Catalog Number: K151KCD-Series
Lot Number: K00E0338
Expiration Date: 30 SEP 2019

Statement: This product was manufactured and tested according to approved MSD documents. This lot meets MSD’s specifications”.

Kit Contents:

Plate					
Lot	Z00E0263	Catalog #	N451KCA-1	Expiration Date:	30 SEP 2019

Antibody Blend	
LotNumber:	D00E0316
Catalog Number:	D21KC-2 / D21KC-3
Expiration Date:	30 SEP 2019
Storage Temperature:	2-8°C
Spot #	Description
1	Human Ang-2

Calibrator Blend	
Lot Number:	A00E0170
Catalog Number:	C01KC-2
Expiration Date:	30 SEP 2019
Storage Temperature	≤ -70°C
Spot #	Description
1	Human Ang-2

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>	08 JUN 2017