



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: MSD® Human IL-6R Ultra-Sensitive Kit

Catalog Number: K151ALC - Series

Lot Number: K0035689

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 #	Z0046772	Catalog #	L451ALA	Expiration Date:	31 AUG 2020

Antibody Blend	
Lot Number:	D0034021
Catalog Number:	D21AL
Expiration Date:	31 AUG 2021
Storage Temperature:	2-8°C
Spot #	Description
1	Anti-hIL-6R Antibody

Calibrator Blend	
Lot Number:	A0032389
Catalog Number:	C01AL
Expiration Date:	31 AUG 2020
Storage Temperature:	≤ -70°C
Spot #	Description
1	Human IL-6R Calibrator

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>	23 APR 2018