



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: MSD® Phospho-VEGFR-2 Whole Cell Lysate Kit

Catalog Number: K151DJD - Series

Lot Number: K0014112

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 #	Z0055920	Catalog #	N451DJB	Expiration Date:	30 NOV 2019

Antibody Blend	
Lot Number:	D0013724
Catalog Number:	D21DJ
Expiration Date:	30 NOV 2019
Storage Temperature:	2-8°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	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	18 OCT 2017