



## CERTIFICATE OF ANALYSIS

COA - 07265

**Product Description:** Phospho-VEGFR-2 Whole Cell Lysate Kit

**Catalog #/Size:** K151DJD-1 / 1 Plate; K151DJD-2 / 5 Plates; K151DJD-3 / 20 Plates

**Lot Number:** K0013793

**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 Number: Z0055403

Plate Lot Catalog Number: N451DJB

Plate Lot Expiration Date: 31 JAN 2018

Antibody blend Lot Number: D0013380

Antibody blend Catalog Number: D21DJ

Antibody blend Expiration Date: 31 AUG 2017

### Specifications:

Release method: Coating uniformity

Mean Intraplate CV:  $\leq 15\%$

Mean Interplate CV:  $\leq 18\%$

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.**

	<b>Name</b>	<b>Function</b>	<b>Signature</b>	<b>Date</b>
<b>Preparer</b>	Xiomara Talavera	Quality Control	<i>X Talavera</i>	02 SEP 2016
<b>Review/Approval</b>	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	02 SEP 2016