



CERTIFICATE OF ANALYSIS

COA - 07723

Product Description: MSD® Human Vascular Injury I Kit

Catalog #/Size: K15135C - Series

Lot Number: K0050731

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: Z0055568

Plate Lot Catalog Number: N45135B

Plate Lot Expiration Date: 31 AUG 2018

Antibody blend Lot Number: D0050666

Antibody blend Catalog Number: D2135

Antibody blend Expiration Date: 31 AUG 2018

Calibrator blend Lot Number: A0020188

Calibrator blend Catalog Number: C0135

Calibrator blend Expiration Date: 31 AUG 2018

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.

	Name	Function	Signature	Date
Preparer	Xiomara Talavera	Quality Control	<i>Talavera</i>	10 JAN 2017
Review/Approval	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	11 JAN 2017