



CERTIFICATE OF ANALYSIS

COA - 07996

Product Description: MSD[®] Human TRAIL Kit

Catalog #/Size: K151KXD-Series

Lot Number: K00E0119

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: Z00E0110

Plate Lot Catalog Number: N451KXA

Plate Lot Expiration Date: 30 JUN 2017

Antibody blend Lot Number: D00E0112

Antibody blend Catalog Number: D21KX

Antibody blend Expiration Date: 30 JUN 2017

Calibrator blend Lot Number: A00E0065

Calibrator blend Catalog Number: C01KX

Calibrator blend Expiration Date: 30 JUN 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.

	Name	Function	Signature	Date
Preparer	Xiomara Talavera	Quality Control		14 MAR 2017
Review/Approval	Vanessa Castleberry	Quality Assurance		15 MAR 2017