



CERTIFICATE OF ANALYSIS

COA - 06294

Product Description: MSD[®] Human Angiopoietin-1 Kit

Catalog #/Size: K151LPD-1 / 1 Plate; K151LPD-2 / 5 Plates; K151LPD-4 / 25 Plates

Lot Number: K00E0258

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: Z00E0207

Plate Lot Catalog Number: N451LPA

Plate Lot Expiration Date: 30 JUN 2018

Antibody blend Lot Number: D00E0232

Antibody blend Catalog Number: D21LP

Antibody blend Expiration Date: 31 MAY 2018

Calibrator blend Lot Number: A00E0122

Calibrator blend Catalog Number: C01LP

Calibrator blend Expiration Date: 30 APR 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	Rolanda Spratley	Biologics/Chemistry		24 MAR 2016
Review/Approval	Dawn Stollar	Quality Assurance		24 MAR 2016