



CERTIFICATE OF ANALYSIS

COA - 06451

Product Description: MSD[®] Human Growth Factor I Kit

Catalog #/Size: K15029C-1/ 1 Plate; K15029C-2 / 5 Plates; K15029C-3 / 20 Plates

Lot Number: K0050620

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: Z0054725
- Plate Lot Catalog Number: N45029B
- Plate Lot Expiration Date: 31 JAN 2017
- Antibody blend Lot Number: D0050547
- Antibody blend Catalog Number: D2029
- Antibody blend Expiration Date: 31 JAN 2017
- Calibrator blend Lot Number: A0020135
- Calibrator blend Catalog Number: C0029
- Calibrator blend Expiration Date: 29 FEB 2016

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		03 FEB 2016
Review/Approval	Dawn Stollar	Quality Assurance		04 FEB 2016