



CERTIFICATE OF ANALYSIS

COA - 06759

Product Description: MSD® Human Osteopontin Kit

Catalog #/Size: K151HJC-1 / 1 Plate; K151HJC-2 / 5 Plates; K151HJC-3 / 20 Plates

Lot Number: K0050713

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: Z0055400
- Plate Lot Catalog Number: N75147A
- Plate Lot Expiration Date: 31 JAN 2018
- Antibody blend Lot Number: D0050639
- Antibody blend Catalog Number: D21HJ
- Antibody blend Expiration Date: 31 JAN 2018
- Calibrator blend Lot Number: A0020177
- Calibrator blend Catalog Number: C0147
- Calibrator blend Expiration Date: 31 JAN 2018

Specifications:

- Release method: Coating uniformity
- Mean Intraplate CV: ≤ 15%
- Mean Interplate CV: ≤ 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		07 APR 2016
Review/Approval	Dawn Stollar	Quality Assurance		07 APR 2016