



## CERTIFICATE OF ANALYSIS

COA - 06752

**Product Description:** MSD<sup>®</sup> Human Alkaline Phosphatase Kit

**Catalog #/Size:** K151HEC-1 / 1 Plate; K151HEC-2 / 5 Plates; K151HEC-3 / 20 Plates

**Lot Number:** K0050581

**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: Z0054439

Plate Lot Catalog Number: N75146A

Plate Lot Expiration Date: 30 APR 2014

Antibody blend Lot Number: D0050502

Antibody blend Catalog Number: D21HE

Antibody blend Expiration Date: 31 JUL 2014

Calibrator blend Lot Number: A0050138

Calibrator blend Catalog Number: C0146

Calibrator blend Expiration Date: 15 Months From Date of Receipt

### 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		07 APR 2016
Review/Approval	Dawn Stollar	Quality Assurance		07 APR 2016