



## CERTIFICATE OF ANALYSIS

COA - 07527

**Product Description:** MSD® Human Myeloperoxidase (MPO) Kit

**Catalog #/Size:** K151EEC-1 / 1 Plate; K151EEC-2 / 5 Plates; K151EEC-3 / 20 Plates

**Lot Number:** K0050723

**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: Z0055486

Plate Lot Catalog Number: L451EEA

Plate Lot Expiration Date: 31 MAR 2018

Antibody blend Lot Number: D0050653

Antibody blend Catalog Number: D21EE

Antibody blend Expiration Date: 31 MAR 2018

Calibrator blend Lot Number: A0020181

Calibrator blend Catalog Number: C01EE

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
<b>Preparer</b>	Xiomara Talavera	Quality Control	<i>X Talavera</i>	23 NOV 2016
<b>Review/Approval</b>	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	28 NOV 2016