



CERTIFICATE OF ANALYSIS

COA - 04238

Product Description: MSD[®] Human Myeloperoxidase (MPO) Kit

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

Lot Number: K0050626

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: Z0054779

Plate Lot Catalog Number: L451EEA

Plate Lot Expiration Date: 31 MAR 2016

Antibody blend Lot Number: D0050556

Antibody blend Catalog Number: D21EE

Antibody blend Expiration Date: 30 APR 2016

Calibrator blend Lot Number: A0020096

Calibrator blend Catalog Number: C01EE

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		18 MAR 2014
Review/Approval	Dana Short	Quality Control		18 APR 2014