



CERTIFICATE OF ANALYSIS

COA - 05667

Product Description: MSD[®] sAPPalpha/sAPPbeta Kit

Catalog #/Size: K15120E-1 / 1 Plate; K15120E-2 / 5 Plates; K15120E-3 / 20 Plates

Lot Number: K0013119

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: Z0054266

Plate Lot Catalog Number: N45120B

Plate Lot Expiration Date: 30 NOV 2013

Antibody blend Lot Number: D0012663

Antibody blend Catalog Number: D21EA

Antibody blend Expiration Date: 30 NOV 2013

Calibrator blend Lot Number: A0010273

Calibrator blend Catalog Number: C01BS

Calibrator blend Expiration Date: 15 Months From Date of Receipt

Calibrator blend Lot Number: A0010295

Calibrator blend Catalog Number: C01BT

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		30 JUN 2015
Review/Approval	Michael Grossman	Quality Assurance		30 JUN 2015