



CERTIFICATE OF ANALYSIS

COA - 06554

Product Description: MSD® sAPPalpha/sAPPbeta Kit

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

Lot Number: K0013812

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: Z0055344
- Plate Lot Catalog Number: N45120B
- Plate Lot Expiration Date: 31 OCT 2017
- Antibody blend Lot Number: D0013384
- Antibody blend Catalog Number: D21EA
- Antibody blend Expiration Date: 31 AUG 2017
- Calibrator blend Lot Number: A0010362
- Calibrator blend Catalog Number: C01BS
- Calibrator blend Expiration Date: 31 JAN 2017
- Calibrator blend Lot Number: A0010363
- Calibrator blend Catalog Number: C01BT
- Calibrator blend Expiration Date: 31 JAN 2017

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	<i>Rolanda Spratley</i>	29 FEB 2016
Review/Approval	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	29 FEB 2016