



CERTIFICATE OF ANALYSIS

COA - 07578

Product Description: MSD® sAPPalpha/sAPPbeta Kit

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

Lot Number: K0013973

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: Z0055683
- Plate Lot Catalog Number: N45120B
- Plate Lot Expiration Date: 31 JAN 2019
- Antibody blend Lot Number: D0013584
- Antibody blend Catalog Number: D21EA
- Antibody blend Expiration Date: 31 JAN 2019
- Calibrator blend Lot Number: A0010391
- Calibrator blend Catalog Number: C01BS
- Calibrator blend Expiration Date: 31 OCT 2018
- Calibrator blend Lot Number: A0010398
- Calibrator blend Catalog Number: C01BT
- Calibrator blend Expiration Date: 31 JAN 2019

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	Xiomara Talavera	Quality Control	<i>Talavera</i>	05 DEC 2016
Review/Approval	Vanessa Castleberry	Quality Assurance	<i>Vanessa Castleberry</i>	06 DEC 2016