



### CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** MSD<sup>®</sup> Human Total PYY Kit

**Catalog Number:** K151MPD - Series

**Lot Number:** K00E0332

**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 #	Z0021262	Catalog #	L15SA	Expiration Date:	30 JUN 2019

Antibody Blend	
LotNumber:	D00E0307
Catalog Number:	N/A
Expiration Date:	31 AUG 2019
Storage Temperature:	2-8°C

Calibrator Blend	
Lot Number:	A00E0165
Catalog Number:	C01MP
Expiration Date:	31 AUG 2019
Storage Temperature:	≤ -70°C

Antibody Blend	
LotNumber:	D00E0308
Catalog Number:	D21MP
Expiration Date:	31 AUG 2019
Storage Temperature:	2-8°C

Specifications:

Release method:

Coating Uniformity	Specifications
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
Review/Approval	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	11 MAY 2017