



## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** MSD® Human FSH Kit

**Catalog Number:** K151ESC - Series

**Lot Number:** K0050796

**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 #	Z0056104	Catalog #	N451ESA	Expiration Date:	30 APR 2020

Antibody Blend	
Lot Number:	D0050737
Catalog Number:	D21ES
Expiration Date:	30 APR 2020
Storage Temperature:	2-8°C

Calibrator Blend	
Lot Number:	A0020231
Catalog Number:	C01ES
Expiration Date:	30 APR 2020
Storage Temperature:	≤ -70°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>	13 MAR 2018