



## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** MSD<sup>®</sup> Human FABP3 Kit

**Catalog Number:** K151HTD-Series

**Lot Number:** K00E0244

**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 #	Z00E0196	Catalog #	N451HTA	Expiration Date:	31 MAR 2018

Antibody Blend	
LotNumber:	D00E0225
Catalog Number:	D21HT
Expiration Date:	30 APR 2018
Storage Temperature:	2-8°C

Calibrator Blend	
Lot Number:	A00E0119
Catalog Number:	C01HT
Expiration Date:	31 MAR 2018
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	Vanessa Castleberry	Quality Assurance	<i>Vanessa Castleberry</i>	18MAY 2017