



### CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** MSD® Human IL-1β Tissue Culture Kit

**Catalog Number:** K151AGB-Series

**Lot Number:** K0035418

**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					
<b>Lot #</b>	Z0046432	<b>Catalog #</b>	L451AGB	<b>Expiration Date:</b>	31 AUG 2019

Antibody Blend	
<b>Lot Number:</b>	D0033819
<b>Catalog Number:</b>	D21AG
<b>Expiration Date:</b>	31 AUG 2019
<b>Storage Temperature:</b>	2-8°C
Spot #	Description
1	Human IL-1beta

Calibrator Blend	
<b>Lot Number:</b>	A0032258
<b>Catalog Number:</b>	C0105
<b>Expiration Date:</b>	31 JAN 2019
<b>Storage Temperature:</b>	≤ -70°C
Spot #	Description
1	Human Tissue Culture Calibrator 2

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
<b>Review/Approval</b>	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	01 FEB 2018