

## CERTIFICATE OF ANALYSIS

Page 1 of 2

**Product Description:** Acute Phase Protein Panel 1 (rat) Kit  
**Kit Catalog Numbers:** K15175C-series  
**Kit Lot Number:** K0040441  
**Expiration Date:** 30 Sep 2020

### Acceptance Data/ Percent Recovery Specifications:

Spot	Name	Calibrator		Control 1		Control 2		Control 3	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	AGP	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
4	A2M	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	AGP	0.146	80% - 120%	< 20%	150	75% - 125%	< 25%
4	A2M	1.00	80% - 120%	< 20%	400	80% - 120%	< 20%

### Acceptance Data/ Non Specific Binding Specifications:

Spot	Name	NSB Spec
1	AGP	< 1.0%
4	A2M	< 1.0%

### Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	AGP	20X Calibrator	8760	ng/mL
		Control 1	54.8	ng/mL
		Control 2	13.7	ng/mL
		Control 3	3.42	ng/mL
4	A2M	20X Calibrator	5540	ng/mL
		Control 1	34.6	ng/mL
		Control 2	8.65	ng/mL
		Control 3	2.17	ng/mL

Control data was generated utilizing internal reference controls.

## CERTIFICATE OF ANALYSIS

Page 2 of 2

Concentrations provided in this table refer to the stock concentration of the calibrators. For top of the standard curve concentrations, please refer to the Standard Curve Data

### Standard Curve Data:


AGP - Spot 1			A2M - Spot 4		
Conc. ng/mL	ECL Signal	%CV	Conc. ng/mL	ECL Signal	%CV
438	91,400	0.9	277	209,000	2.9
110	63,200	3.1	69.3	71,200	4.0
27.4	30,300	4.3	17.3	17,600	2.4
6.84	9,720	3.1	4.33	4,300	3.4
1.71	2,750	2.5	1.08	1,310	2.0
0.428	893	2.5	0.271	581	2.4
0.107	415	3.9	0.0676	399	2.3
0.00	258	6.0	0.00	349	2.4

### Components:

Description	Lot Number	Storage Temperature	Expiration Date
Acute Phase Protein Panel 1 (rat) Plate	Z0056307	2-8°C	31 Jan 2021
Anti-rat AGP Antibody	D0040801	2-8°C	30 Apr 2022
Anti-rat A2M Antibody	D0040800	2-8°C	31 Jan 2022
Acute Phase Protein Panel 1 (rat) Calibrator Blend	A0040219	≤ -70°C	31 Mar 2021
Diluent 100	Y0270255	2-8°C	30 Sep 2020
MSD Read Buffer T	Not Kit Specific	Not Kit Specific	N/A

### Statement:

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		05 DEC 2018