

## CERTIFICATE OF ANALYSIS

Page 1 of 2

**Product Description:** Acute Phase Protein Panel 1 (rat) Kit  
**Kit Catalog Numbers:** K15175C-Series  
**Kit Lot Number:** K0040461  
**Expiration Date:** 31 Mar 2021

**Acceptance Data/ Percent Recovery Specifications:**

Spot	Name	Calibrator		Acute Phase Control 1		Acute Phase Control 2	
		Recovery	CV	Recovery	CV	Recovery	CV
1	AGP	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
4	A2M	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
		Acute Phase Control 1	1010	ng/mL
		Acute Phase Control 2	89.2	ng/mL
4	A2M	20X Calibrator	5540	ng/mL
		Acute Phase Control 1	569	ng/mL
		Acute Phase Control 2	51.5	ng/mL

Control data was generated utilizing internal controls.

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

## CERTIFICATE OF ANALYSIS

Page 2 of 2

### Standard Curve Data:


AGP - Spot 1			A2M - Spot 4		
Conc. ng/mL	ECL Signal	%CV	Conc. ng/mL	ECL Signal	%CV
438	177,000	2.2	277	206,000	1.4
110	133,000	2.3	69.3	76,400	2.5
27.4	61,300	1.9	17.3	20,500	2.3
6.84	17,800	1.8	4.33	5,180	3.3
1.71	4,470	2.8	1.08	1,610	2.5
0.428	1,280	2.0	0.271	659	1.9
0.107	477	2.4	0.0676	433	2.3
0.00	213	3.8	0.00	347	2.1

### Components:

Description	Lot Number	Storage Temperature	Expiration Date
Acute Phase Protein Panel 1 Plate	Z0056482	2-8°C	31 Oct 2021
Anti-rat AGP Antibody	D0040830	2-8°C	31 Aug 2022
Anti-rat A2M Antibody	D0040829	2-8°C	31 Aug 2022
Acute Phase Protein Panel 1 (rat) Calibrator Blend	A0040219	≤ -70°C	31 Mar 2021
Diluent 100	Y0270262	2-8°C	30 Sep 2021
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>	Karen Hamilla	Quality		06 AUG 2019