

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: Acute Phase Protein Panel 1 (rat) Kit

Kit Catalog Numbers: K15175C-Series

Kit Lot Number: K0040478

Expiration Date: 31 Mar 2021

Acceptance Data/ Percent Recovery Specifications:

Spot	Name	Calibrator		Acute Phase Protein Panel 1 (rat) Control 1		Acute Phase Protein Panel 1 (rat) Control 2		Acute Phase Protein Panel 1 (rat) 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%	220	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 Protein Panel 1 (rat) Control 1	110	ng/mL
		Acute Phase Protein Panel 1 (rat) Control 2	27.4	ng/mL
		Acute Phase Protein Panel 1 (rat) Control 3	6.84	ng/mL
4	A2M	20X Calibrator	5540	ng/mL
		Acute Phase Protein Panel 1 (rat) Control 1	69.3	ng/mL
		Acute Phase Protein Panel 1 (rat) Control 2	17.3	ng/mL
		Acute Phase Protein Panel 1 (rat) Control 3	4.33	ng/mL

Control data was generated utilizing internal reference 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	163,000	1.4	277	143,000	2.2
110	122,000	2.1	69.3	53,000	2.9
27.4	58,200	3.4	17.3	14,200	3.8
6.84	18,100	3.2	4.33	3,630	5.1
1.71	4,850	3.8	1.08	1,090	3.7
0.428	1,500	2.7	0.271	451	2.3
0.107	622	4.5	0.0676	292	1.6
0.00	348	4.0	0.00	236	4.0

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Acute Phase Protein Panel 1 Plate	Z0056554	2-8°C	31 Dec 2021
Anti-rat AGP Antibody	D0040876	2-8°C	31 Dec 2022
Anti-rat A2M Antibody	D0040875	2-8°C	31 Dec 2022
Acute Phase Protein Panel 1 (rat) Calibrator Blend	A0040219	≤ -70°C	31 Mar 2021
Diluent 100	Not Kit Specific	Not Kit Specific	N/A
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
Review/Approval	Xiomara Talavera	Quality		01 SEP 2020