

CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Acute Phase Protein Panel 1 (rat) Kit
Kit Lot Number: K0040509
Expiration Date: 30 Apr 2023
Kit Catalog Numbers: K15175C-Series

Precision and Accuracy Specifications:

Spot	Name	Calibrator		Acute Phase Control 1		Acute Phase Control 2		Acute Phase 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%	90.6	80% - 120%	< 20%

Non-Specific Binding Specifications:

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

Stock Concentration of Calibrator:


Spot	Name	Conc.	Units
1	AGP	11620	ng/mL
4	A2M	7020	ng/mL

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Acute Phase Protein Panel 1 Plate	Z0056818	2-8°C	30 Apr 2023
Anti-rat AGP Antibody	D0040939	2-8°C	31 May 2024
Anti-rat A2M Antibody	D0040938	2-8°C	31 May 2024
Acute Phase Protein Panel 1 (rat) Calibrator Blend	A0040257	≤ -70°C	30 Jun 2023

Additional components may be provided to execute the assay. Additional components are not lot-specific.

The above product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		05 MAR 2021