



EXTENSION MEMO


Page 1 of 1

Product Description(s): Reference Standard 1

Catalog Number(s) / Size(s): C00ADK-2

Lot Number: A0080286

Based on additional testing, the original date of expiration for this Reference Standard has been extended by 60 months. The expiration date of A0080286 has been extended to 28 FEB 2031.

Function	Name	Signature	Date
Quality	Karen Hamilla		06 MAY 2024

This memo supersedes all other statements regarding date of expiration for this product.



RFE MEMO

Page 1 of 1

Product Description(s): Serology Control Pack 1


Catalog Number(s) / Size(s): C4381-1

Lot Number: A00C0825, A00C0826, A00C0827

Based on additional testing, the original date of expiration for this Control Pack has been extended by 60 months. The expiration date of A00C0825, A00C0826, A00C0827 has been extended to 31 MAR 2032

The expiration dates for the following controls have been extended:

Description	Lot Number	Original Expiration Date	New Expiration Date
Serology Control 1.1	A00C0825	31 MAR 2027	31 MAR 2032
Serology Control 1.2	A00C0826	31 MAR 2027	31 MAR 2032
Serology Control 1.3	A00C0827	31 MAR 2027	31 MAR 2032

Function	Name	Signature	Date
Quality	Karen Hamilla		03 MAY 2024

This memo supersedes all other statements regarding date of expiration for this product.



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: V-PLEX Respiratory Panel 6 (IgG) Kit, SECTOR
Lot Number: K0082831
Expiration Date: 31 MAR 2028
Catalog Numbers: K15755D-Series

Kit Components:

The kit expiration date provided in the COA is determined by the component with the earliest expiration date. As the kit is supplied with different catalog products, not all configurations include all the components. In such cases, the applicable expiration date is determined by the shortest expiry among the components included in the kit that is received.

Description	Lot Number	Storage Temperature	Expiration Date
Respiratory Plate 6 96-Well 10-Spot SECTOR Plate	Z0057615	2-8°C	31 Mar 2028
SULFO-TAG MSD GOLD Anti-Human IgG Antibody	D00V0043	2-8°C	31 Jul 2028
Reference Standard 1	A0080286	≤ -70°C	28 Feb 2031
Serology Control 1.1	A00C0825	≤ -70°C	31 Mar 2032
Serology Control 1.2	A00C0826	≤ -70°C	31 Mar 2032
Serology Control 1.3	A00C0827	≤ -70°C	31 Mar 2032
Diluent 100	Not Kit Specific	2-8°C	See Label
MSD Blocker A Kit	Not Kit Specific	Room Temperature	See Label
MSD GOLD Read Buffer B	Not Kit Specific	Room Temperature	See Label

Functional Testing Results:

Spot	Sample Type	Calibrator	Controls		Samples	
	Metric	Avg Signal Ratio Test:Reference	Control Recovery	Control Conc. CV	Avg % Difference Test:Reference	Slope Test:Reference
1	Flu A/Wisconsin/2019 H1	50 – 200%	70 – 130%	< 20%	± 30%	0.70 – 1.3
2	Flu A/Darwin/2021 H3	50 – 200%	70 – 130%	< 20%	± 30%	0.70 – 1.3
3	Flu B/Austria/2021 HA	50 – 200%	70 – 130%	< 20%	± 30%	0.70 – 1.3
6	hPIV1 F	50 – 200%	70 – 130%	< 20%	± 30%	0.70 – 1.3
7	hPIV2 F	50 – 200%	70 – 130%	< 20%	± 30%	0.70 – 1.3
8	hPIV3 F	50 – 200%	70 – 130%	< 20%	± 30%	0.70 – 1.3
9	hPIV4 F	50 – 200%	70 – 130%	< 20%	± 30%	0.70 – 1.3

Additional Comments:

Functional testing was executed using Respiratory Plate 6 (Z0057615), SULFO-TAG Anti-Human IgG Antibody (D00V0043), Reference Standard 1 (A0080286), and Serology Controls (A00C0825, A00C0826, A00C0827).

All kit components were manufactured and tested according to MSD documents. The lots listed in the Kit Components table meet MSD's specifications and plate spot layout.

Statement:

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		18 NOV 2025