

# AcroMetrix™ EDTA Plasma Dilution Matrix 45 mL

For research use only. Not for use in diagnostic procedures.

**REF** S2284 AcroMetrix EDTA Plasma Dilution Matrix 45 mL

## Product Use

AcroMetrix™ EDTA Plasma Dilution Matrix 45 mL, a human plasma matrix shown to be non-reactive for blood borne pathogens, viral nucleic acids and serological markers, is intended for use in preparing specimens for testing in plasma based nucleic acid assays. AcroMetrix EDTA Plasma Dilution Matrix is not intended for use as a substitute for the integral diluents, standards or controls provided by in vitro diagnostic kit manufacturers. For Research Use Only. Not for use in diagnostic procedures.

## Summary and Explanation

AcroMetrix EDTA Plasma Dilution Matrix 45 mL is an EDTA based normal human plasma (NHP) that has been tested and found non-reactive for HBV DNA, HCV RNA, HIV-1 RNA, anti-HIV-1/2, HBsAg, anti-HBc, anti-HCV and anti-HTLV I-II.

AcroMetrix EDTA Plasma Dilution Matrix 45 mL provides manufacturing firms, clinical laboratories and blood banks with well-characterised, pre-screened negative diluent. The use of a standard dilution matrix also allows testing laboratories and other end users to compare results obtained at different times, by different operators, or by different methodologies, and ensures that the sample diluent itself is not contributing to testing variability.

## Principles of the Procedure

AcroMetrix EDTA Plasma Dilution Matrix 45 mL has been carefully formulated to mimic naturally occurring normal human plasma. Therefore, the product can be used with any test procedure that requires preliminary dilutions of test specimens in human serum or plasma.

## Reagents

Catalog Number	Product Name	Quantity
S2284	AcroMetrix EDTA Plasma Dilution Matrix 45 mL	1 x 45 mL

## Precautions and Warning

**WARNING:** Although the NHP used in the production of this product was determined to be non-reactive for antibodies to HIV-1 and HIV-2, HIV-1 RNA, HBsAg, antibodies to HCV, HCV-RNA, HBsAg, antibodies to HBc, HBV-DNA, and antibodies to HTLV I-II, it should be handled as if capable of transmitting infectious agents. Contains ≤0.05% Sodium azide.

Observe universal precautions for prevention of transmission of infectious agents when handling these materials<sup>1,2,3</sup>.

Do not pipette by mouth. Use personal protective equipment, including lab coats, gloves and safety glasses. Do not eat, drink or smoke in areas where diluents and specimens are handled. Disinfect liquids, materials or spills with a 0.5% sodium hypochlorite solution. Dispose of all materials and liquids used in the procedure as if they contained pathogenic agents.

## Storage Instructions

It is recommended that AcroMetrix EDTA Plasma Dilution Matrix 45 mL be stored at -20°C or lower to ensure highest quality. AcroMetrix EDTA Plasma Dilution Matrix may be refrozen after first use, thawed and used a second time without risk to product quality. Discard any unused material after the second use.

## Instructions for Use

Thaw AcroMetrix EDTA Plasma Dilution Matrix 45 mL quickly at 37°C to reduce the risk of cryoprecipitate formation, vortex briefly, and store at 2 to 8°C until ready to use. Return any unused matrix to the recommended storage conditions immediately after use.

AcroMetrix EDTA Plasma Dilution Matrix 45 mL should be handled and tested in a manner identical to that required for serum/plasma diluents or clinical specimens utilized in the nucleic acid test procedure being evaluated. Follow the manufacturer's instructions and recommendations for the handling and testing of diluents and clinical specimens.

## Limitations

AcroMetrix EDTA Plasma Dilution Matrix 45 mL is intended for use as a diluent for use in routine laboratory testing procedures where patient specimens may require dilution. It is not intended for use as a substitute for the integral diluents, standards or controls provided by *in vitro* diagnostic kit manufacturers. This product is for Research Use Only. Not for use in diagnostic procedures.

## References

- Centers for Disease Control (CDC). Recommendations for prevention of HIV transmission in health care settings. MMWR 1987; 36 (supplement no. 2S).
- Centers for Disease Control (CDC). Update: Universal precautions for prevention of transmission of human immunodeficiency virus, hepatitis B virus, and other bloodborne pathogens in health-care settings. MMWR 1988; 37:377-388.
- Centers for Disease Control (CDC). Guidelines for prevention of transmission of human immunodeficiency virus and hepatitis B virus to health-care and public-safety workers. MMWR 1989; 38(S-6): 1-36.

## Glossary:

<http://www.thermofisher.com/symbols-glossary>



Microgenics Corporation  
46500 Kato Road  
Fremont, CA 94538 USA  
US Customer and  
Technical Support:  
1-800-232-3342



For insert updates go to:  
[www.thermofisher.com/Acrometrix](http://www.thermofisher.com/Acrometrix)