

AcroMetrix™ PeliSpy™ Sero Chagas Positive Control

thermo
scientific

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

REF 956805 AcroMetrix PeliSpy Sero Chagas Positive Control

Product Use

The AcroMetrix™ PeliSpy™ Sero Chagas Positive Control is used to compare precision and accuracy of results between laboratories and among lots of commercial assays and for the identification of variability with the diagnostic method. For Research Use Only. Not for use in diagnostic procedures.

Composition

The AcroMetrix PeliSpy Sero Chagas Positive Control contains a pool of high titer anti-*Trypanosoma cruzi* sera in a normal human serum (NHS) matrix. The diluent matrix is comprised of defibrinated, delipidized normal human plasma with preservatives, tested and found negative for HIV-1 RNA, anti-HIV-1, anti-HIV-2, anti-HTLV I – II, HBV DNA, HBsAg, HCV RNA, anti-HCV, anti-HBs, and anti-*T. cruzi*.

Principles of the Procedure

The AcroMetrix PeliSpy Sero Chagas Positive Control has been carefully formulated to mimic naturally occurring human specimens containing anti-*T. cruzi*. Therefore, the controls can be used with any test procedure designed for detecting anti-*T. cruzi* in human serum or plasma.

Expected results utilizing the AcroMetrix PeliSpy Sero Chagas Positive Control must be established by the end user for their anti-*T. cruzi* assay.

Reagents

Catalog Number	Control Name	Quantity
956805	AcroMetrix PeliSpy Sero Chagas Positive Control	10 x 3.5 mL

Precautions and Warning

WARNING: The AcroMetrix PeliSpy Sero Chagas Positive Control contains anti-*T. cruzi* and should be considered potentially biohazardous. Observe universal precautions for prevention of transmission of infectious agents when handling this material^{1, 2, 3}.

H412 - Harmful to aquatic life with long lasting effects.

Avoid release to the environment. Dispose of contents/container to location in accordance with local/regional/national/international regulations.

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 controls 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.

This product contains ProClin™ 300 as a preservative. Use caution when working with and disposing of this material.

Storage Instructions

It is recommended that the AcroMetrix PeliSpy Sero Chagas Positive Control be stored at 2°C to 8°C to ensure highest quality. Once opened, the vial of AcroMetrix PeliSpy Sero Chagas Positive Control can be used up to four (4) weeks from date of opening if stored according to instructions.

Instructions for Use

Bring the AcroMetrix PeliSpy Sero Chagas Positive Control to room temperature just before use and mix tubes by gently swirling. Return to proper storage conditions immediately after use.

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

© 2018 Thermo Fisher Scientific Inc. All rights reserved. Proclin is a trademark of Rohm and Haas Company. All other trademarks are the property of Thermo Fisher Scientific or its affiliates.

MAN0004347-5
2018 06