## Not for Clinical Use.

REF 956980 AcroMetrix BCR-ABL Panel Kit

AcroMetrix™ BCR-ABL Panel is intended for being used as an external control panel for analytical evaluation of BCR-ABL test methods. The product is not intended for clinical use.

### **Summary and Explanation**

AcroMetrix BCR-ABL Panel includes 5 panel members. The product contains lyophilized human cell lines K-562 (BCR-ABL e14a2 fusion gene positive) cells and HL-60 (BCR-ABL negative) cells mixed at different ratios. Each vial contains approximately 1 X 10<sup>6</sup> cells.

AcroMetrix BCR-ABL Panel is traceable to international standard (IS) and is manufactured using similar approaches to the 1<sup>st</sup> World Health Organization International Genetic Reference Panel for Quantitation of BCR-ABL Translocation by RQ-PCR (WHO Primary Reference Panel NIBSC 09/138) (White et al., 2010)

### **Contents and Traceability**

Catalog Number	Product Name		Quantity
956980	BCR-ABL Panel		1 vial each x 5
Panel Member		Targeted Value <sup>IS</sup>	
Panel A		10% BCR-ABL/ABL	
Panel B		1% BCR-ABL/ABL	
Panel C		0.1% BCR-ABL/ABL	
Panel D		0.01% BCR-ABL/ABL	
Panel E		0.0032% BCR-ABL/ABL	

### Warnings and Precautions

AcroMetrix BCR-ABL panel contains biological material of human origin. As with all the samples of biological origin, AcroMetrix BCR-ABL Panel should be handled and discarded according to guidelines issued by your institution's health, safety and environment office.

AcroMetrix BCR-ABL Panel is contained in glass vials. Care should be taken during handling to avoid breakage and injuries inflicted upon by broken glass.

### Instructions for Preparation

- 1. RNA extraction should be performed using your local standard procedures for the extraction of RNA from human blood samples.
- Do not attempt to remove any portion of the lyophilized material from the vials prior to reconstitution.
- 3. Vials are single use only. Extract RNA from the entire content of each vial.
- 4. Vials used in the same study on the same date should be processed simultaneously.
- 5. Lysate of the lyophilized material should be processed immediately.
  - It is preferable to inject by RNase-free syringe the required volume of cell lysis buffer prior to opening the vials, to avoid loss of material due to pressure inside the vial.
  - Use a volume appropriate for 1 x 10<sup>6</sup> cells. Add the lysis buffer directly onto the lyophilized material inside the vial.
  - Unscrew the cap and set aside. Keep the rubber stopper on.
  - Using a syringe fitted with a needle, aspirate the desired volume of lysis buffer from its container.
  - · Inject the lysis buffer through the rubber stopper.
  - Replace the cap and mix by gently inverting a few times till the lysis buffer is clear and no particulate material is visible.
  - Before opening, tap the vials gently to collect the lysed materials at the bottom of the vial.
  - Thoroughly mix by pipetting the lysis buffer up and down for at least 10 times.
  - Process the entire lysate immediately according to your local RNA extraction procedures.

### **Storage Condition**

All vials should be stored immediately at -20°C upon receipt.

### Reference

- WHO International Standard 1<sup>st</sup> WHO International Genetic Reference Panel for quantitation of BCR-ABL translocation by RQ-PCR (NIBCS code: 09/138), Instructions for use, Version 4.0.
- 2. White H.E. et al, Establishment of the 1st World Health Organization International Genetic Reference Panel for quantitation of BCR-ABL mRNA. Blood, 2010; 116(22): 3111-7.

### **Glossary:**

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

# 444

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