# PanoHealth® Dried Blood Sample Elution Buffer

For eluting a dried blood sample

Catalog #: PANO-DBSEB

Instructions for Use Last Revised Jun. 10<sup>th</sup>, 2020

Caution: Extraordinary useful information enclosed



ISO 13485:2016

3607 Parkway Lane, Suite 100 Peachtree Corners, GA 30092

Tel: 1-888-494-8555 (Toll Free) or 770-729-2992, Fax:770-206-2393 Web: www.RayBiotech.com, Email: info@raybiotech.com

## I. INTRODUCTION

Dried blood samples are routinely used for the collection, testing, and archiving of blood samples (capillary blood, venous blood, serum, or plasma). Collected samples can be analyzed for proteins, antibodies, DNA, or other biomarkers. The PanoHealth® Dried Blood Sample Elution Buffer is used for optimal elution of the sample for downstream analysis using a number of applications. This product is for research use only. Please read the instructions carefully before starting your experiment

## II. MATERIAL PROVIDED

- 1. 1X Phosphate Buffered Saline with Tween® 20 6 mL
- 2. Protease Inhibitor Cocktail 1 vial of lyophilized powder

# III. ADDITIONAL MATERIAL REQUIRED

- 1. Precision pipettes to deliver 20 µl to 1 ml volumes
- 2. Orbital Shaker
- 3. Microcentrifuge tubes
- 4. Microcentrifuge
- 5. Vortex

#### IV. STORAGE

Upon receipt, the 1X Phosphate Buffered Saline with Tween® 20 should be stored at room temperature. The Protease Inhibitor Cocktail should be stored at -80°C until reconstitution immediately before use.

## V. ELUTION BUFFER PREPARATION

- 1. Briefly spin down the Protease Inhibitor Cocktail vial
- 2. Add 100 µL of 1X Phosphate Buffered Saline with Tween® 20 into the Protease Inhibitor Cocktail Vial and pipette up and down to mix thoroughly.
- 3. Transfer the  $100 \,\mu\text{L}$  of reconstituted Protease Inhibitor Cocktail from step 2 back into the 1X Phosphate Buffered Saline with Tween® 20 bottle. Mix well.

# VI. DRIED BLOOD SAMPLE ELUTION PROCEDURE

- 1. Elute the dried blood sample at a 1:10 ratio. For example: 1 5x10mm strip of the PanoHealth® Blood Collection Device (<u>PANO-BC-CD</u>) should be eluted with 250 μL of Elution Buffer.
- 2. Elute for 4 hours at room temperature on an orbital shaker, vortexing for 10 seconds every 30 minutes
- 3. Transfer the supernatant to a clean tube leaving the filter paper behind
- 4. Centrifuge at 14,000 rpm for 10 minutes
- 5. Collect the supernatant into a clean tube
- 6. Following elution, the eluate should be stored at 4°C for no more than 24 hours; eluates should be aliquoted, labeled, and stored at -80°C until use.

This product is for research use only.



©2020 RayBiotech, Inc.