



EAGLE
BIOSCIENCES

Calprotectin Sample Collection Kit

Catalog Number: CAL35-C50

For Research Use Only. Not for use in diagnostic procedures.

v. 12 (30 SEP 23)

EAGLE BIOSCIENCES, INC.
20A Northwest Blvd., Suite 112, Nashua, NH 03063
Phone: 617-419-2019 Fax: 617-419-1110
WWW.EAGLEBIO.COM



INTENDED USE

The Eagle Biosciences Calprotectin Sample Collection Kit is a device intended for human fecal sample collection and extraction. The extracted sample can be used directly for calprotectin level determination by professional laboratories. The Calprotectin Sample Collection Kit is for research use only and not to be used for diagnostic procedures.

For further information about this kit, its application, or the procedures in this insert, please contact the Technical Service Team at Eagle Biosciences, Inc at www.EagleBio.com or at 866-411-8023.

MATERIALS PROVIDED

1. Fecal Calprotectin Sample Collection Tube: 48 units/kit. Each tube contains 4.5 mL Calprotectin fecal sample collection and extraction buffer.
2. Instructions for Use

SAFETY PRECAUTIONS

The reagents must be used in a professional laboratory environment and are for research use only. Wear gloves when performing sample collection and treat the reagents as if they are hazardous. The buffer may cause irritation on skin upon contact. Do not get in the eyes, on skin, or clothing. Never transfer any reagents with mouth. Do not ingest or inhale. On contact, flush with copious amounts of water for at least 15 minutes. Exercise Good Laboratory Practices.

MATERIALS REQUIRED BUT NOT PROVIDED

1. Plastic wide-mouth stool collection container

SPECIMEN COLLECTION AND PREPARATION

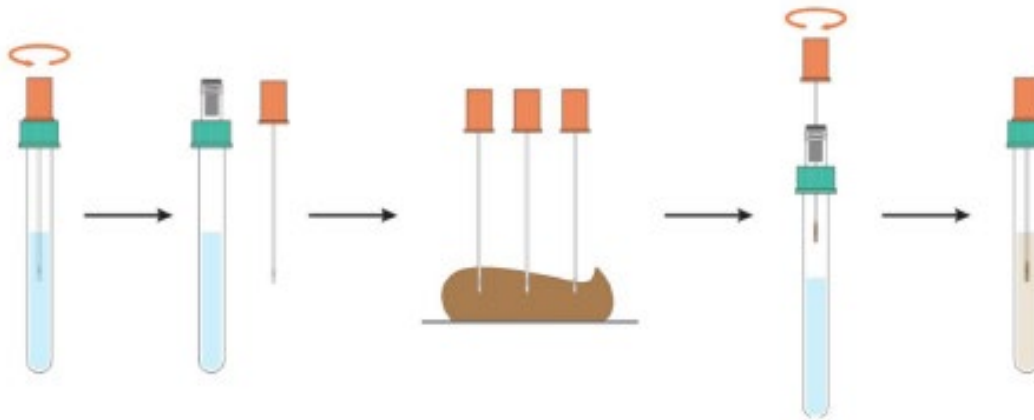
1. Open the lid of primary stool collection container.
2. Open the device sampling lid (orange color) and pull the sampling lid out of the tube.
3. Use the grooved region of the wand to stir and mix soft stool or dip into hard stool multiple times (3 – 5 sites) to collect a small amount of stool sample.
4. Insert the sampling wand and screw the sampling lid back into the sampling tube in a vertical position. Tightly seal the sampling lid against the tube. Let the sample soak in extraction buffer at least 10 minutes.
5. Vortex the tube on upright position at least 10 seconds to dissolve the stool sample completely. (a digital roller mixer can be used to assist in facilitating the mixing process)
6. The collected sample should be on upright position on a rack for 1 – 5 minutes to sediment stool particles before assay.

PROCEDURAL NOTES

1. This device is recommended for the secondary collection of stool sample in professional laboratories, while patient should perform primary collection of stool sample into a plastic wide-mouth stool collection container in bathroom.
2. Careful technique is necessary to ensure reproducibility and maximum detective sensitivity.
3. For liquid stool please use pipette to add 100µL sample directly into Fecal Calprotectin Sample Collection Tube.



SAMPLE COLLECTION PROCEUDURE



STORAGE

The extracted stool sample should be tested withing 8 hours, or stored at 2-8oC and tested within 3 days. It may be stored below -20oC for a longer storage period. Avoid more than three freeze – thaw cycles for each specimen.

WARRANTY INFORMATION

Eagle Biosciences, Inc. warrants its Product(s) to operate or perform substantially in conformance with its specifications, as set forth in the accompanying package insert. This warranty is expressly limited to the refund of the price of any defective Product or the replacement of any defective Product with new Product. This warranty applies only when the Buyer gives written notice to the Eagle Biosciences within the expiration period of the Product(s) by the Buyer. In addition, Eagle Biosciences has no obligation to replace Product(s) as result of a) Buyer negligence, fault, or misuse, b) improper use, c) improper storage and handling, d) intentional damage, or e) event of force majeure, acts of God, or accident.

Eagle Biosciences makes no warranties, either expressed or implied, except as provided herein, including without limitation thereof, warranties as to marketability, merchantability, fitness for a particular purpose or use, or non-infringement of any intellectual property rights. In no event shall the company be liable for any indirect, incidental, or consequential damages of any nature, or losses or expenses resulting from any defective product or the use of any product. Product(s) may not be resold, modified, or altered for resale without prior written approval from Eagle Biosciences, Inc.

For further information about this kit, its application or the procedures in this kit, please contact the Technical Service Team at Eagle Biosciences, Inc. at info@eaglebio.com or at 866-411-8023.