



DCM063-3
Ed. 03/2012

SALIVA COLLECTION DEVICE

IVD

LOT

See external label

2°C 8°C



Σ = 100 PZ

REF DKO063

INTENDED USE

Eagle Biosciences [Saliva Collection Device](#) is for collection of saliva samples using the glass collection device.

1.1 Reagent and material supplied in the kit

1. Glass Collection Tubes 100 pz **REF DCE063/6301-0**
2. Plastic Straw 100 pz **REF DCE0063/6302-0**

2.1 Method

Collect saliva samples at the times indicated.

If no specific instructions have been given, the following should be noted:

- a) **If saliva is collected in the morning this should be carried out before you brush your teeth.**
- b) **During the day do not have anything to eat or drink for at least one hour before you take your saliva sample**
- c) **To ensure the best quality of samples, it is advisable that samples are collected and sent to the Laboratory as early in the week as possible – this will avoid any delays in a post over a weekend.**
- d) **It is very important that a good clear sample is received – i.e. no contamination from food, lipstick, blood (bleeding gums) or other matter.**

3.1 Sample collection instructions

- 1) Carefully remove the red top from the glass tube.
- 2) Using the straw provided, gently expel saliva through the straw and into a glass tube.
- 3) Ensure that the tube is **at least half full of saliva (NB: bubbles do not count!)** – this may take any time up to 10 minutes – **the amount and the quality of the sample are both important.**
- 4) If you have difficulty activating saliva, you may chew gently a chewing gum.
- 5) Remove the straw and discard.
- 6) Firmly replace the top on each glass tube.

3.2 Technician instructions

- 1) **Once received the saliva tube (with sample) centrifuge for 15 minutes at 3000 rpm.**
- 2) **Store at – 20°C for at least 1 hour.**
- 3) **Centrifuge again for 15 minutes at 3000 rpm.**
- 4) **The saliva sample is ready to be tested.**
- 5) **Store the sample at 2- 8°C for one week or at – 20°C for longer times.**

Ed. 03/2012

DCM063-3

DiaMetra S.r.l. Headquarter: Via Garibaldi, 18 – 20090 SEGRATE (MI) Italy

Tel. 0039-02-2139184 – 02-26921595

Fax 0039-02-2133354.

Manufactory: Via Pozzuolo 14, 06038 SPELLO (PG) Italy

Tel. 0039-0742-24851

Fax 0039-0742-316197

E-mail: info@diametra.com

distributed in the US/Canada by:

EAGLE BIOSCIENCES, INC.

20A NW Blvd, Suite 112 Nashua, NH 03063

Phone: 617-419-2019 FAX: 617-419-1110

www.EagleBio.com

• info@eaglebio.com



EAGLE
BIOSCIENCES