



**G-Biosciences ♦ 1-800-628-7730 ♦ 1-314-991-6034 ♦ [technical@GBiosciences.com](mailto:technical@GBiosciences.com)**

A Geno Technology, Inc. (USA) brand name

# SalivaOptimize™ Proteinase K

For saliva based SARS CoV-2 testing  
(Cat. # 786-1783, 786-1784)



**think proteins! think G-Biosciences [www.GBiosciences.com](http://www.GBiosciences.com)**

## INTRODUCTION

SalivaOptimize™ Proteinase K has been optimized to provide maximum proteolytic activity within a short period for rapid and maximum release of Covid-19 viral nucleic acid for RT-PCR detection technique.

## ITEM(S) SUPPLIED

Cat. #	Description	Size
786-1783	SalivaOptimize™ Proteinase K [20mg/ml]	1 ml
786-1784	SalivaOptimize™ Proteinase K [20mg/ml]	10 ml

## STORAGE CONDITION

The product is shipped with blue ice. Upon arrival store at -20°C

## SPECIFICATIONS

- **Form:** liquid
- **Activity:**  $\geq 30$  U/mg
- **Specific Activity:**  $\geq 40$  U/mg protein
- DNase and RNase free (Molecular grade)

## PREPARATION BEFORE USE

Use SalivaOptimize™ Proteinase K [20mg/ml] directly as indicated in protocol. Make single use aliquots, store frozen and avoid repeated thawing.

## PROTOCOL

**Saliva collection Precautions:** Saliva should be collected with the assistance of trained healthcare worker or technician. For self-collection, get approval from appointed diagnostic lab for Covid-19 testing

1. Clean the hands with sanitizer or soap and water. Wear appropriate PPE.
2. Ensure that collection container is properly labeled with sample identifying information.
3. Direct the sample provider (patient) to begin pooling saliva in mouth.  
**NOTE:** Only natural saliva which naturally pools in mouth is collected. No coughing or/and sniffing prior to sample collection is required.  
**NOTE:** Water must be avoided 10 minutes before sample collection. Other foods and drinks should be avoided 30 minutes before sample collection.
4. Remove the lid of the container and ask the patient/sample provider to gently pour the saliva into the container until the volume reaches 0.5ml.  
**NOTE:** Volume exclude bubbles
5. Close the lid and clean the outer surface of the container with 70% ethanol or 10 % bleach.
6. Proceed to sample processing.

### **Saliva storage/transportation**

Saliva sample can be processed immediately. If transportation to other facility is required, it can be shipped at room temperature as virus in saliva is stable room temperature for at least 7 days. Once sample reaches in laboratory it can be processed immediately or store at 4°C for short term (up to 7 days from day collection) or -80°C for long term storage.

### **Proteinase K treatment of saliva**

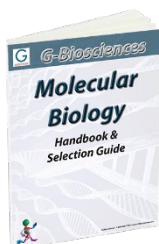
1. Assemble the required number of tubes (PCR or other types) on rack or on the automated platform.
2. Aliquot 5 µl SalivaOptimize™ Proteinase K per tube.
3. Vortex the saliva samples till they are homogenous.
4. Transfer 50µl of saliva sample per tube containing proteinase K solution.
5. Close the tubes lids tightly and vortex for 1 minute at 3000-5000 rpm.
6. Briefly centrifuge the tubes.
7. Incubate the samples for 5 minutes at room temperature
8. Heat the samples at 95°C for 5 minutes on PCR machine or appropriate thermocycler.
9. Briefly centrifuge the tubes to collect sample at the bottom of the tube.
10. Store the processed sample at -80°C or use immediately for RT-qPCR testing.
11. Use 5 µl of this Proteinase K treated saliva sample for RT-qPCR.

### **STORAGE**

Store SalivaOptimize™ Proteinase K solution in small aliquots at -20°C or -80°C.

### **RELATED PRODUCTS**

Download our Molecular Biology Handbook.



<http://info2.gbiosciences.com/complete-molecular-biology-handbook>

For other related products, visit our website at [www.GBiosciences.com](http://www.GBiosciences.com) or contact us.



**[www.GBiosciences.com](http://www.GBiosciences.com)**