

Diagenode One

Small, Light, Perfect

Diagenode Sonication Machine



Dear customer,

Thank you for choosing our device and for your confidence in Diagenode.

We have designed this device using our highest level of knowledge in shearing so that it exceeds your expectations. This device will serve your research needs for many years and will provide the best experience with NGS.

INTENDED USE

The Diagenode One is intended for laboratory use. It is not suitable for clinical use. Use of the Diagenode One in any manner other than as directed herein could cause harm to persons and may void the warranty. Diagenode will not be responsible for injury or damage resulting from improper use of the Diagenode One.

WARNING

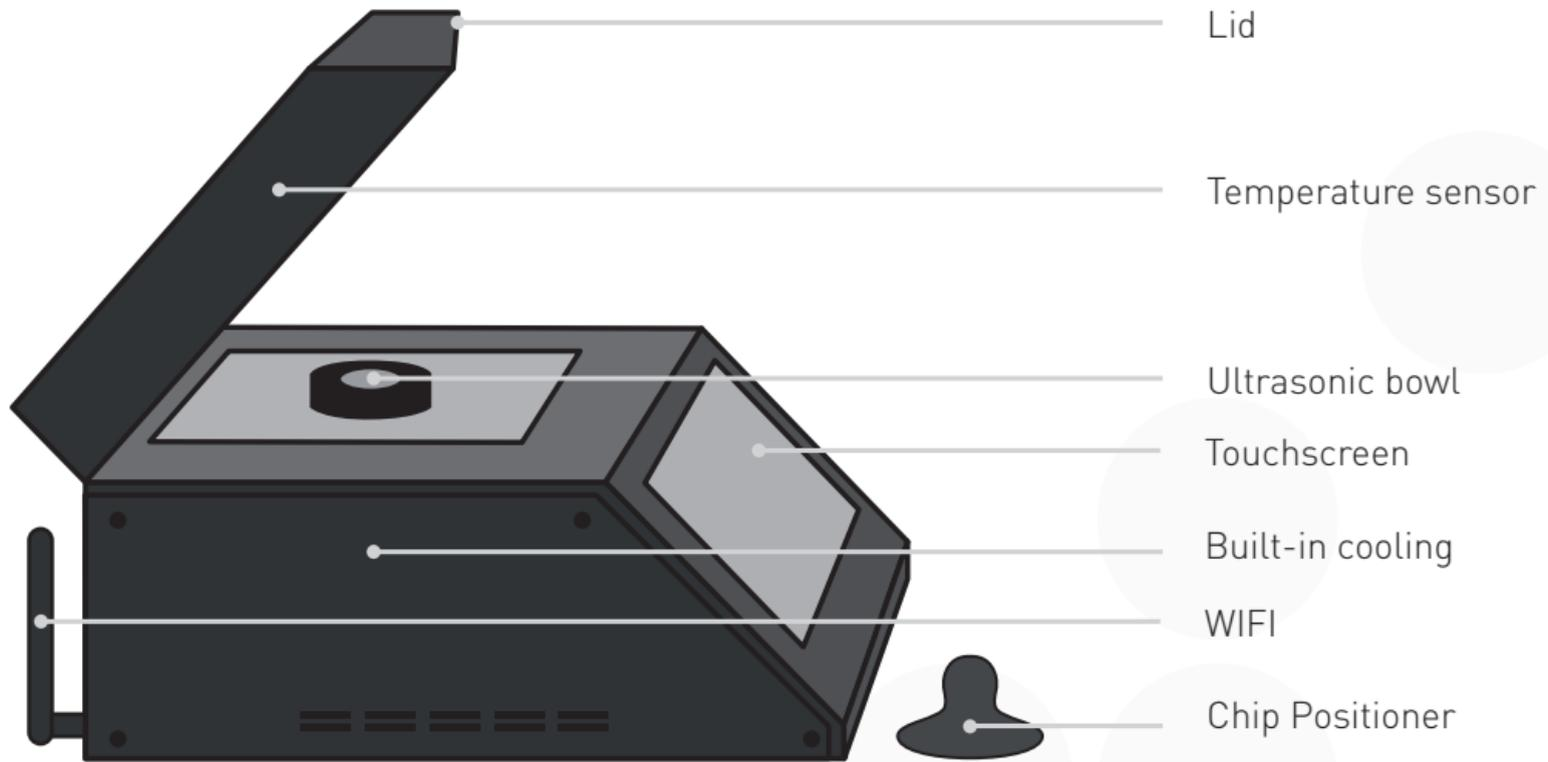
Carefully read the following instructions before using your Diagenode One to avoid the risk of personal injury and damage.

- Do not use the Diagenode One for anything other than its intended use.
- After removing the packaging, make sure that the Diagenode One is intact.
- Do not allow the Diagenode One to be used by untrained individuals.
- Do not start the sonication if the gel and the chip are not in position.
- Clean the Diagenode One carefully and regularly with soft tissue and water.
- Do not use any accessories not recommended by Diagenode.
- Do not use the Diagenode One if it is showing any sign of breakdown or malfunction.
- Do not attempt to dismantle or repair.
- Keep far grid clear on the back panel.

Description

TECHNICAL SPECIFICATIONS

Input voltage	100-240
Maximum consumed current	1.56 A
Unit dimension	200 (W) x 285 (D) x 165 (H) mm
Total weight	5 kg
Disposable	Micro fluidic chip + ultrasonic gel



Description

INTERFACE



Sonication : Running a shearing experiment (follow instructions)



Time : Display the current time and date



Temperature : Display the temperature at the ultrasonic bowl



Settings : Visualize and change the parameters of the Diagenode One



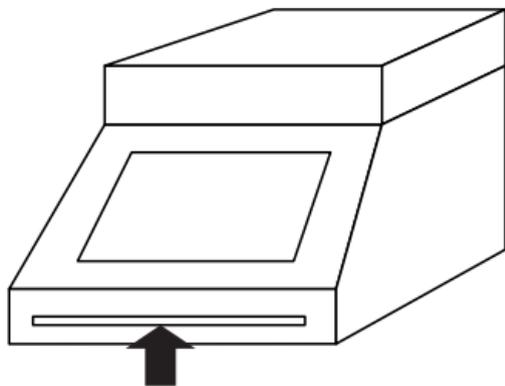
Info : General information about your Diagenode One



Manual : Operation instructions for best DNA shearing with the Diagenode One

LED LIGHT BAR

For best results, the Diagenode One is equipped with a fully integrated cooling system. The perfect control of the temperature provides optimal DNA shearing and sample quality protection. The Diagenode One operates through sonication and cooling cycles. The LED light bar allows the tracking of the processing.



The **blue light** indicating cooling time.

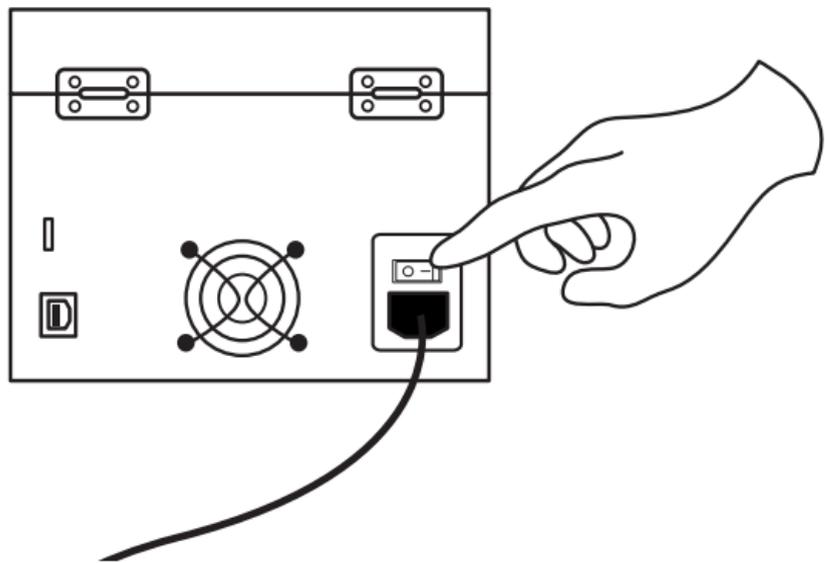
The **red light** corresponding to sonication.

The **green light** indicates the end of the shearing.

In addition, the LED lights indicate the progression time.

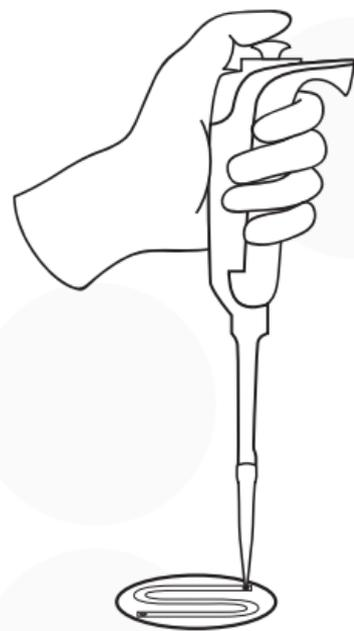
1

Start the Diagenode One and let it cool down
(optional change the set temperature)



2

Load the microfluidic chip with your sample



Instructions

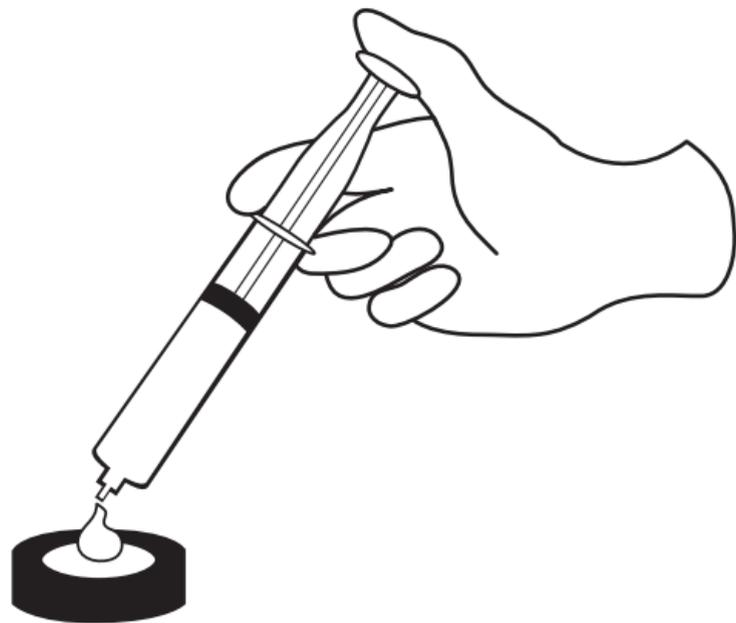
3

Cover the microfluidic chip with the tape and **press** to ensure a good bond



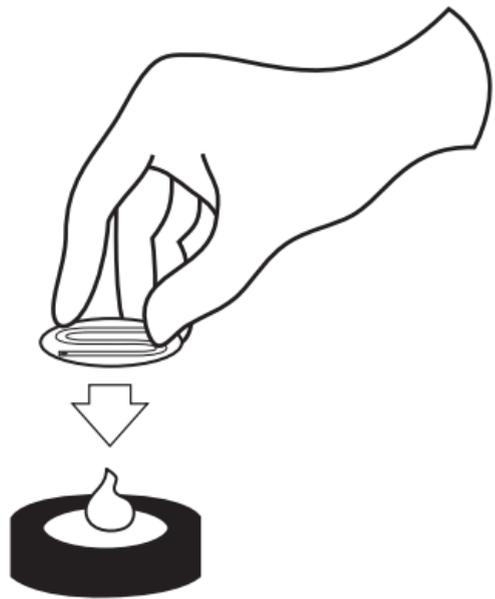
4

Dispense 1ml of gel on the ultrasonic bowl



5

Install the chip (**holes face up**) off at the ultrasonic bowl



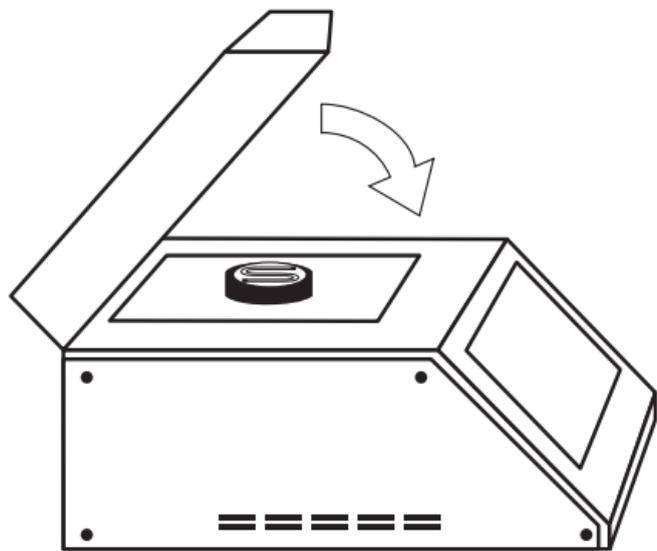
6

Gently **press** the chip with the **chip positioner** to set a good position

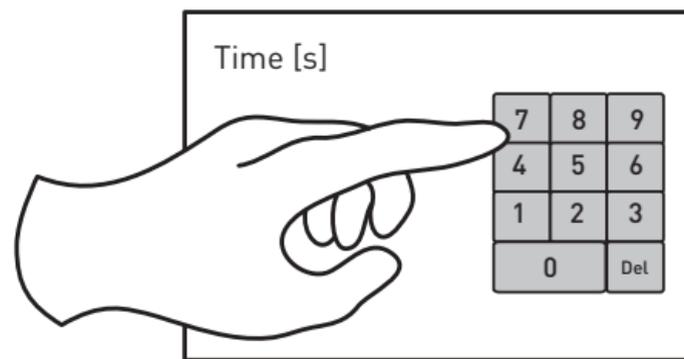


Instructions

7 Close the lid

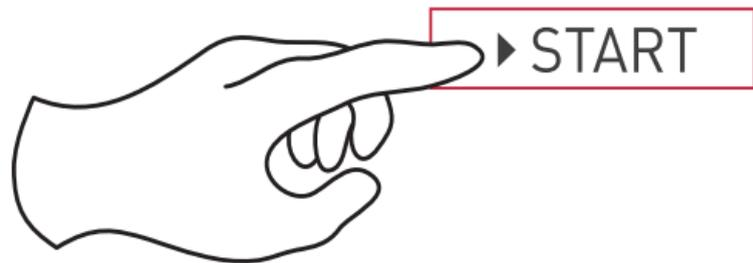


8 Set sonication time*



*Visit <https://www.diagenode.com/en/dna-shearing-guide> to find the adequate length of sonication for your expected DNA size

9 Start the sonication



10 Collect your sample and clean the ultrasonic bowl with a soft cloth or with a paper towel



CLEANING

Clean the sonication area to remove all remaining gel. For all others areas, we highly recommend using only a humid soft cloth when cleaning.

SAFETY APPROVAL

Diagenode One has been assessed and **certified under the CE standards.**

ENVIRONMENTAL PROTECTION

Packaging materials contain recyclable materials and the Diagenode One has been manufactured using high quality materials that can be recycled.

WARRANTY

Diagenode guarantees all products from any manufacturing defects as we rigorously test all products to meet strict quality standards. Diagenode warrants that all standard components of its instruments will be free of defects in materials and workmanship for a period of one (1) year from the date that the warranty period begins, unless the original quotation or accompanying documentation states a different warranty period. All warranty periods begin on the date of delivery and apply only to the first purchaser of the product. If a manufacturing defect arises and a valid claim is received within the warranty period, Diagenode, at its discretion, will repair or replace the product in accordance with the warranty terms and conditions stated herein. In case of repair or replacement of a product under warranty, Diagenode will cover the expenses to return the repaired or replacement product.

This warranty covers only manufacturing defects and does not cover any damage caused by misuse, lack of compliance to recommendations stated in the manual, neglect, accidents, abrasion, or exposure to extreme temperatures, chemical solvents, or acids. Improper or incorrectly performed maintenance or repairs will void the warranty.



© 2016 Diagenode, Inc. All rights reserved. The content of this document cannot be reproduced without prior permission of the authors.