

# **Thermo Scientific Nalgene Magnetic Carboy Stirrer**

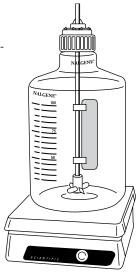
#### **General Information**

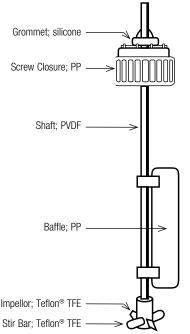
Thermo Scientific™ Nalgene™ Magnetic Carboy Stirrer is a magnetic stir bar suspended from a shaft for use with a magnetic stir plate. This reusable accessory is designed to stir the contents of a 10-L or 20-L carboy with 83B (83mm) neck finish.

**Note:** This accessory is not designed for cell culture use.

Nalgene<sup>™</sup> Magnetic Carboy Stirrer, Cat. No. DS2227-0020, consists of:

- 83B (83mm) screw closure, polypropylene
- Grommet, silicone
- Shaft, PVDF with caps (stainess steel 304-reinforced interior)
- Baffle, polypropylene
- Stir bar, Teflon® TFE
- Impeller, Teflon® TFE





## Assembly

This product is supplied assembled. Two additional impellers have been supplied as replacement parts.

#### Use

The Nalgene Magnetic Carboy Stirrer may be adjusted for use with either a 10-L or 20-L carboy. to adjust the magnetic carboy stirrer height, simply push the shaft through the grommet in the top of the closure to the approximate height of the carboy. When the magnetic carboy

stirrer is installed in the carboy, the stir bar should sit approximately 1" 925.4mm) from the bottom of the carboy.

Insert the Magnetic Carboy Stirrer into the carboy through the carboy neck opening. You may flex the impeller ends together for an easier fit through the opening.

The stir speed depends upon the magnetic stir plate used, the viscosity of the solution being stirred, and the distance between the stir bar and the magnetic stir plate. Maximum stirring efficiency can be achieved in two ways:

- 1. Adjust the stir control on the magnetic stir plate
- 2. Adjust the baffle on the magnetic carboy stirrer: the baffle prevents whirlpool and vortex formation. Move the baffle along the shaft's length for optimum mixing.

#### **Chemical Resistance**

The Nalgene Magnetic Carboy Stirrer is resistant to strong acids, stong bases, and alcolhols.

## Cleaning

Use a warm solution of a mild, neutral detergent. We recommend Thermo Scientific Nalgene L900 liquid detergent (Cat. No. 900-4000). Rinse thoroughly with distilled water. Air dry if possible.

#### **Autoclaving**

The Nalgene Magnetic Carboy Stirrer can be autoclaved within a carboy.

#### Autoclaving the stirrer within a carboy

#### **CAUTION**

Before autoclaving, make sure that the carboy is autoclavable. Check the resin code on the bottom of the carboy. DO NOT autoclave HDPE or LDPE carboys.

- Clean the carboy and magnetic stirrer and rinse thoroughly with distilled water, being sure to rinse off any trace chemicals or detergent.
- 2. Adjust the height of the stirrer so it does not touch the bottom of the carboy. Place the magnetir carboy stirrer inside the carboy.
- Before autoclaving, make sure that the threads are disengaged.
   Magnetic stirrer closures must be completely disengaged from the carboy neck threads for proper steam sterilization to take place and to prevent implosion of the carboy.
- 4. Closure and mgnetic stirrer shaft should sit completely vertical when the threads are disengaged.
- 5. Autoclave for 20 minutes at 121°C, 15 psig (1.0 bar).



# **Thermo Scientific Nalgene Magnetic Carboy Stirrer**

#### Autoclaving the stirrer by itself

- Clean the carboy and magnetic stirrer and rinse thoroughly with distilled water, being sure to rinse off any trace chemicals or detergent.
- 2. Autoclave for 20 minutes at 121°C, 15 psig (1.0 bar).

#### DO NOT exceed these recommendations

### **Temperature Limits**

The polypropylene (PP) screw closure and baffle can withstand temperatures to 135°C.

The Teflon® TFE stir bar and impeller can withstand temperatures to  $121^{\circ}\text{C}$ .

The silicone grommet can withstand temperatures from -20°C to 121°C.

The PVDF shaft can withstand temperatures to 121°C.

## Replacement Part

| Part                          | Cat. No.     |
|-------------------------------|--------------|
| 83B Screw Closure and Grommet | 71-2227-1020 |

Visit www.thermoscientific.com product resources pages for complete product use guidelines.

For more information, see the current Lab Consumables catalog, or contact Technical Support at: **technical.nalgene@thermofisher.com** 

