

Instruction Manual Note Vs2 (22.04.2008)

Diagenode sa

CHU, Tour GIGA, 3e étage
Avenue de l'hôpital, 1
4000 Liège - BELGIUM
Tel. +32 (0)4 364 20 50
info@diagenode.com
www.diagenode.com

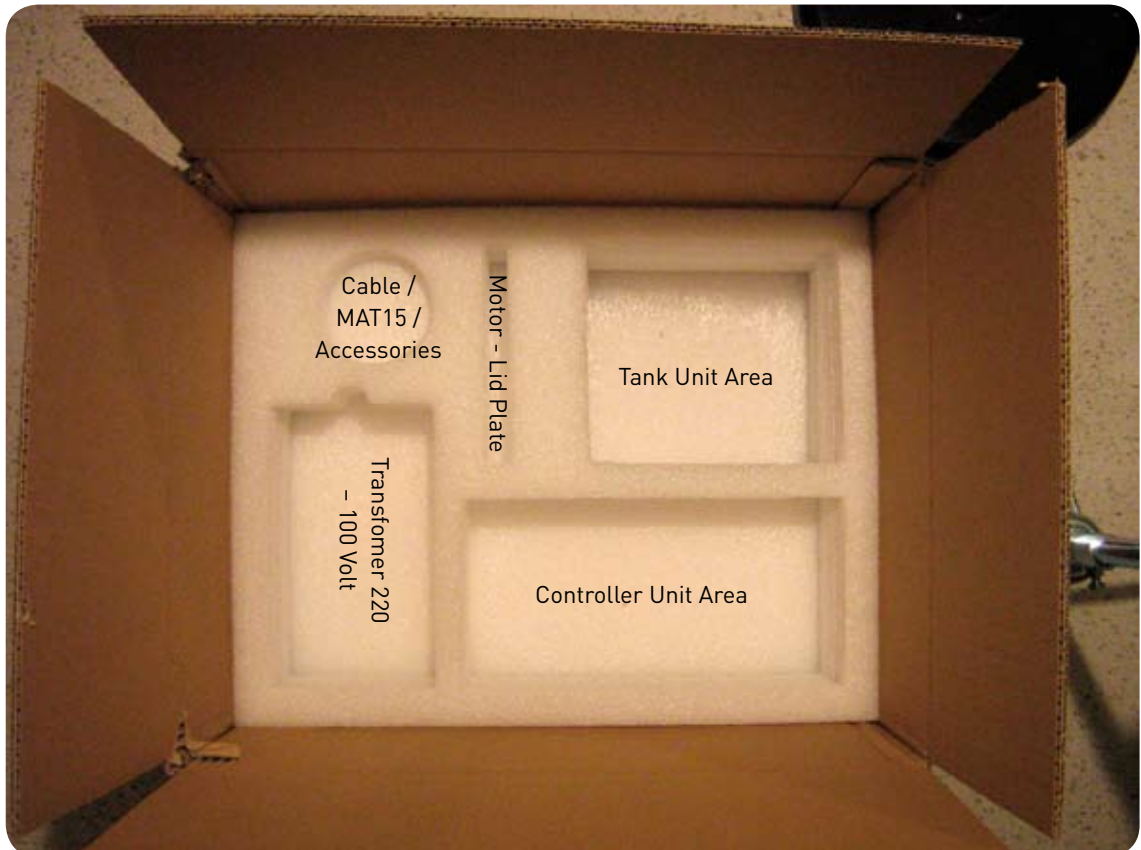
Bioruptor™

User manual

Packaging Instructions

Due to the extreme fragility of Bioruptors it is critical that they be packaged correctly. The following are steps in correct packaging as shown, using our pre-moulded container and their components.

The box is composed of a double wall corrugated carton, a foam base and a foam lid made of ultra shock-absorbent foam. Each part of Bioruptor has a designated area in the box.



1. Tank unit preparation

- Place a scrap piece of corrugated cardboard over the top of the tank (as shown).
- Tie packaging strap around the tank and cardboard.
- Tighten the strap as securely as possible (due to extreme fragility of tank unit which is prone to separating damage during transit).



2. Placement of components

- The controller unit should be placed in the largest rectangular recessed area in the moulded foam box with the fan toward the outer side of the box (as shown).
- The tank unit should be placed in the square recessed area in the moulded foam box towards the inside of the box (as shown). Wedge a piece of cardboard in the space between the housing and the foam to keep the tank in place.
- The motor plate should be placed in the slotted area in the moulded foam, the motor towards the inside of the box (as shown). Tuck the cable into the foam in the middle of the circle of the motor plate to block its movement.
- Roll the grey cable up and place it in the circular recessed area in the moulded foam box. Place the Mat-15 and other accessories in the same circular area.
- The transformer (220V-100V) unit should be placed in the remaining rectangular recessed area in the moulded foam box with the cable between the foam and housing (as shown). Note: The transformer is not used in every area of the world.



3. Placement of foam lid and seal box

- Place foam lid on the top so that the protrusion in the lid is over the transformer's area (as shown). (The protrusion helps to keep things in place during transit.)
- The carton box should be easy to close; otherwise you might have to shift the Bioruptor parts and wires slightly.
- Close the box and seal securely with wide tape.





Instruction Manual (version 01)

Diagenode sa
CHU, Tour GIGA, 3e étage
Avenue de l'hôpital, 1
4000 Liège - BELGIUM
Tel. +32 (0)4 364 20 50
info@diagenode.com
www.diagenode.com