

PRODUCT NAME	
Metallic soundproof box	
Cat. No: <b>B01200001</b> (BioBox-200)	Format: 1 unit

**Product description**

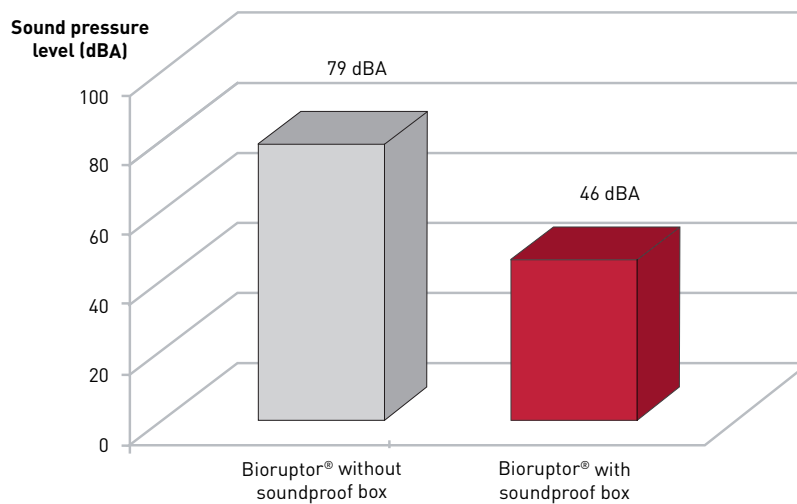
This metallic soundproof box has been designed for a highly efficient acoustic insulation of the Bioruptor® sonication bath. This box is only compatible with Bioruptor® Standard and Bioruptor® Plus. Can be stored at room temperature or in a cold room (4°C).

**Dimensions :**

Width : 35 cm  
Depth : 35 cm  
Height : 52 cm



The metallic soundproof box reduces dramatically the acoustic intensity level of the Bioruptor®.



**Figure 1.** Sound pressure of the Bioruptor® without and with metallic soundproof box has been measured in a soundproof room.