



HyPerforma Single-Use Mixer DS 300

Uncrating Instructions

© 2013 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific Inc. and its subsidiaries. Specifications, terms and pricing are subject to change. Not all products are available in all countries. Please consult your local sales representative for details.

Initial Setup Instructions

The S.U.M. DS 300 will arrive with the following items:

- Docking Station mixer unit
- Hand-held control device
- Motor and mount
- Optional drum positioner
- Drive shafts (customer-specified)
- Torque and spanner wrenches
- Equipment Turnover Package (ETP), the standard or detailed documentation package (if ordered) that is located on the USB drive provided with mixer

S.U.M. DS 300 drums will arrive with the following items:

- Outer support container or drum
- Drum Dolly

Hardware Uncrating

The Single-Use Mixer DS 300 will arrive crated. Be sure to follow the unpacking instructions provided.

IMPORTANT NOTE: When uncrating hardware, retain all packing materials for possible future use.

Follow the steps below to unpack the S.U.M. DS 300.

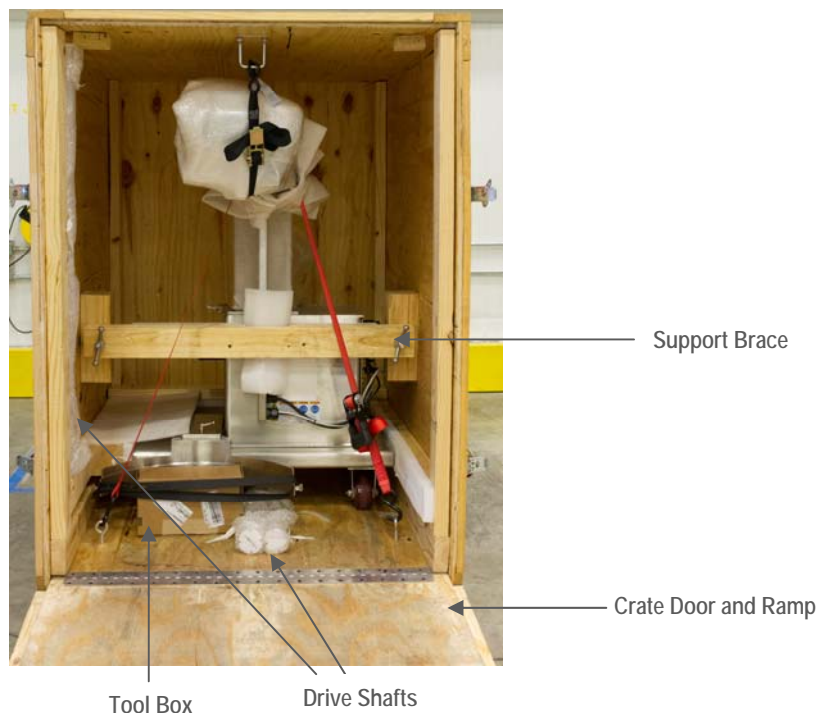
1. Remove the security screws from each of the four latches on either side of the front of the crate as shown in Figure 1, below.

Figure 1: Removing security screws from crate latches



2. Swing the crate door down. The door will create a ramp for moving items in and out, as shown in Figure 2, below.

Figure 2: Open crate



3. Remove the drive shafts from the crate and set them aside. Depending on the size and number of drive shafts ordered, they may be secured to the bottom of the crate. Longer shafts will be secured upright in the front corner of the crate.
4. Remove the uncrating and crating instructions from the clear plastic holder affixed to the inside wall of the crate.
5. Remove the small cardboard box that contains the tools. Unwrap the tools and set them aside.
6. Remove the strap that secures the mixer to the top of the crate, as shown in Figure 3.

Figure 3: Removing the top strap



7. Remove the strap that secures the S.U.M. DS 300 to the eyes screwed into the bottom of the crate, as shown in Figure 4.

Figure4: Removing the bottom strap



8. Remove the large wing nuts from the front support brace as shown in Figure 5. Then pull the front support brace from the crate as shown in Figure 6. Pull both sides straight out at the same time to keep the bolts from binding inside the brace.

Figure 5 and 6: Removing the wing nuts and support brace



9. Remove the S.U.M. DS 300 unit carefully, by rolling it forward out of the crate, as shown in Figure 7.



WARNING: If you need to lift the unit, lift it by the bottom of the cart frame. **Do not lift the unit by its casters.**

Figure 7: Rolling the Docking Station out of the crate



10. Remove all foam and plastic wrap packaging materials from the unit. The hand-held control unit is wrapped separately and attached to the side of the handle. Take care when cutting the packaging material from this device. Then hang the device on the hook as shown in Figure 8.

Figure 8: Hanging the hand held device on its hook



11. If your unit is equipped with the optional rotating motor head, rotate the motor into working position as shown in Figure 9.

Figure 9: Motor head rotation (if unit was ordered with this option)



12. Inspect the unit to ensure that no damage occurred during shipping.

NOTE: Contact your sales representative immediately if damage has occurred.

13. Place all packaging and shipping materials back into the crate for possible future use.
14. Place the bearing hub, torque wrench, and spanner wrench into the appropriate storage holes on the Docking Station unit. See Figure 10.

Figure 10: Stored bearing hub and wrenches



15. Place the drive shaft(s) in the appropriate storage holder on the Docking Station. See Figure 11.

Figure11: Drive shafts stored in the Docking Station



UK/Ireland +44 1442 233555
USA and Canada +847 295 7500

Thermo
S C I E N T I F I C
Part of Thermo Fisher Scientific