

Technical Bulletin

Strip-Fuge™ Operating Instructions

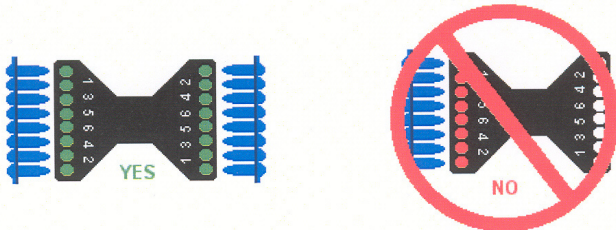
Technical Bulletin AL-254



1. Place the Strip-Fuge on a flat, clear surface.
2. To open, press firmly down on the release tab while lifting the dome.
3. Close dome and activate side-switch. Device will not operate with dome open.
4. To stop, deactivate side-switch. Wait until rotor is completely stopped before releasing the dome. Remove samples.

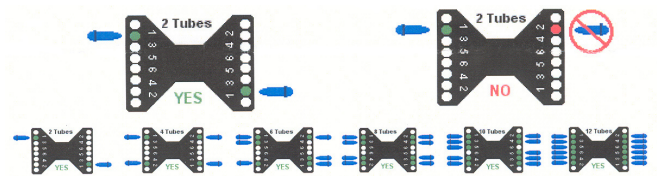
Loading the Butterfly Rotor 16 x 0.2 mL Tube Strips

- Load all tubes on both strips with equal amounts of sample. Close the dome securely.
- Insert strips into rotor. If necessary, add a strip of dummy-tubes to provide balance.

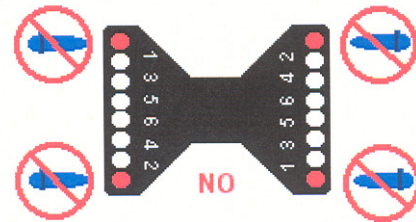


Loading the Butterfly Rotor 0.2 mL Single Tube Strips

- Load tubes with equal amounts of sample in opposing positions as indicated on the rotor.



- Never place single tubes in corner positions. Your unit has been designed to minimize footprint. Single tubes may make contact with the dome and cause damage if placed in corner positions.



Safe Handling

1. Do not immerse device in water. Injury due to shock or fire may result.
2. Do not operate device with dome open, or open dome before rotor has ceased spinning. Do not insert non-standard tubes or other foreign articles into rotor.
3. Do not operate under uneven loads. Destabilized operation may result in damage or injury.
4. Store unit at room temperature in a dry area. Avoid sun, moisture or extreme temperatures.

Sigma brand products are sold through Sigma-Aldrich, Inc. Sigma-Aldrich, Inc. warrants that its products conform to the information contained in this and other Sigma-Aldrich publications. Purchaser must determine the suitability of the product(s) for their particular use. Additional terms and conditions may apply. Please see reverse side of the invoice or packing slip.