

# GeneChip Hybridization Oven 645i: best practices

This quick reference card was developed to aid in the proper use of the Applied Biosystems™ GeneChip™ Hybridization Oven 645i.

## Recommendations

Never stack more than two ovens.



Place ovens away from air conditioning units.

Ensure the proper amount of free space around the oven.



Do not attempt to monitor the thermal performance of the oven using temperature probes other than what is provided in the kit. The temperature profile is calibrated at the factory to facilitate proper operation.

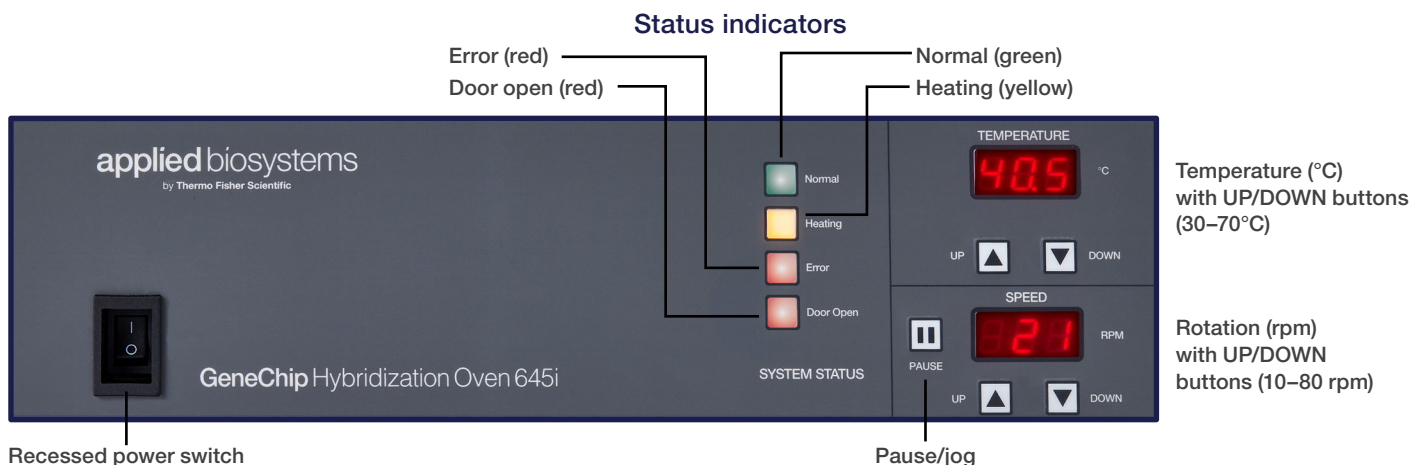





Allow for adequate air flow, i.e., do not block the rear vent.



Balance the carriers in the rotisserie by placing them 180° from each other.

Use the carrier retainer to secure the cartridge carriers.



| If you want to:  | Then do this:   | Description   |
|--|---|---|
| Turn on/off power  | <ul style="list-style-type: none"> <li>Press the I/O button on the front/left.</li> <li>Press <b>PAUSE</b> to view the current settings and operate the oven.</li> </ul>  | <p>This button controls the line power to the unit.</p>           |
| Adjust the oven temperature  | <ul style="list-style-type: none"> <li>Press the <b>UP</b> or <b>DOWN</b> button to set the temperature.</li> </ul>   | <p>These controls are associated with the heating function.</p>  |
| Adjust the rotisserie motor speed, pause the motor, or jog the motor | <ul style="list-style-type: none"> <li>Press and hold the <b>UP</b> or <b>DOWN</b> button until the set point blinks to access the set point.</li> <li>Press the <b>UP</b> or <b>DOWN</b> button to set the speed (rpm).</li> <li>Press <b>PAUSE</b> to pause the rotisserie.</li> <li>When the door is open, you can use the <b>PAUSE</b> button as the <b>JOG</b> button.</li> <li>Press <b>PAUSE</b> again to take the oven out of pause to program and run the oven.</li> </ul> | <p>These controls adjust the speed of the rotisserie motor.</p>  |