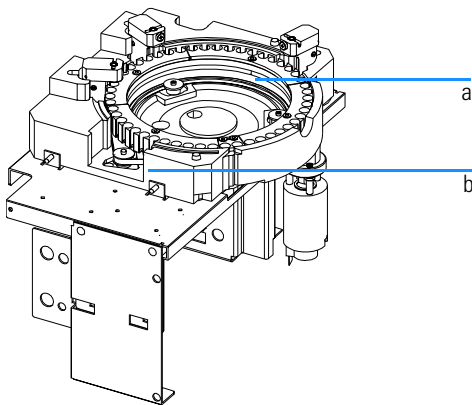




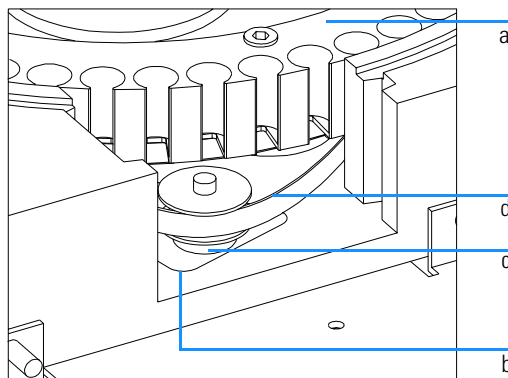
Installing the Tray Belt Modification (G1600-60026)

- 1** Remove tray base assembly.
Remove foam parts (a) and foam insert (b)
from the sample carousel.

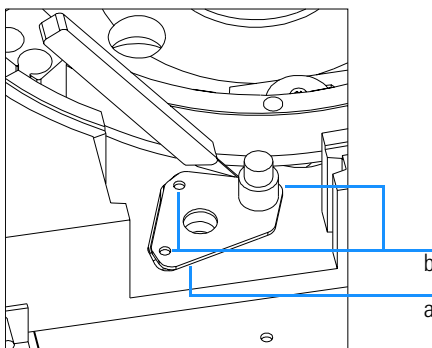
Do not remove the heat exchanger.



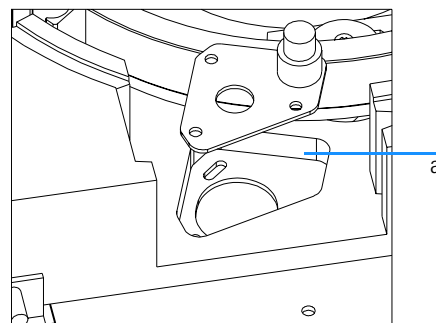
- 2** Remove all segments (a).
Loosen screws that hold tray motor (b).
Remove motor tooth wheel by opening the
two small hex screws (c).
Remove tray motor and tray belt (d).



- 3** Put roller plate on top of foam (a).
Make sure the three holes fit correctly (b).
Use a knife to cut away overlapping foam.

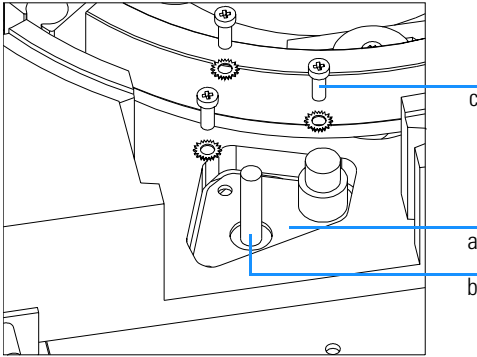


- 4** The new cut-out must look like this (a).
Remove all foam cuttings. Remaining cuttings
inhibit proper function.

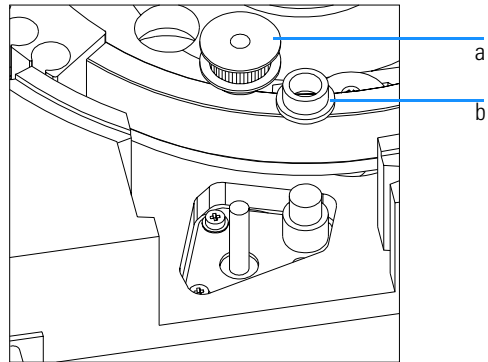


Installing the Tray Belt Modification (G1600-60026)

- 5** Insert the roller plate (a).
Put the tray motor back (b).
Use the provided screws and fix the motor (c). Do not tighten it completely, because the tension of the tray belt needs to be adjusted



- 6** Put the motor tooth wheel (a) back and tighten the two hex screws.
Insert the bearing sleeve (b).

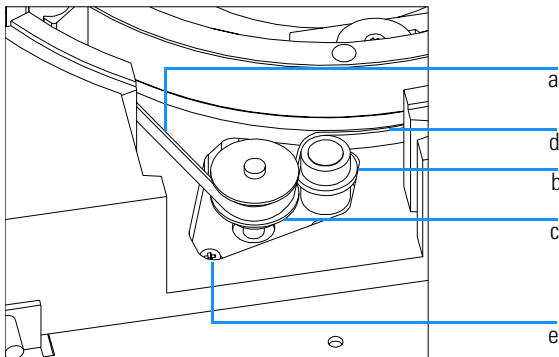


- 7** Install the longer tray belt (a) part number 1500-1040, labeled B340 MXL.

Ensure proper alignment of the tray belt on the belt roller (b), the tooth wheel (c) and the tray (d).

Move the motor to give the tray belt a certain tension. It is not necessary to over do it.

Tighten the tray motor screws (e).



G1600-90109

© Copyright Agilent Technologies 2000

Printed in Germany Edition 02/00

Part No. G1600-90109