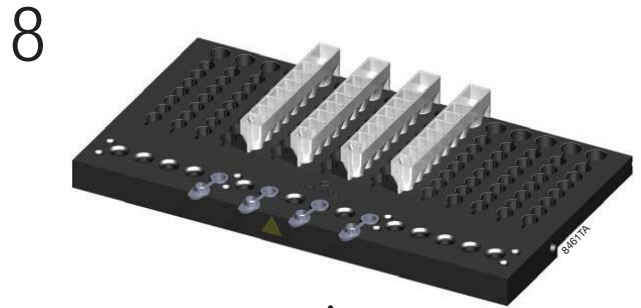
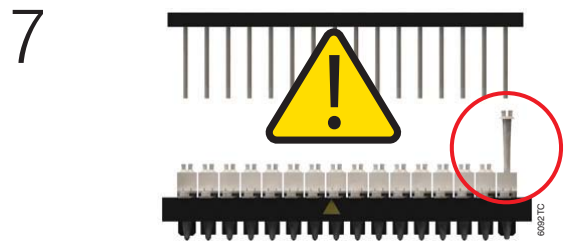
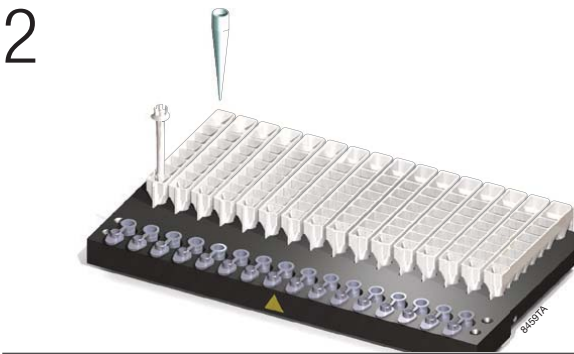




i Please see Maxwell 16 Technical Manual for detailed instructions.

Instructions for Use of Maxwell[®] 16 Instrument.
Configured in LEV.



Cleaning



Turn off and unplug instrument.



1. Extend platform and wipe off with 70% ethanol then dH₂O.



2. Wipe down magnetic rods and plunger bar adaptor with 70% ethanol.



3. Wipe underside of plunger bar with 70% ethanol then dH₂O.



4. Clean platform slides with a dry cloth.



Promega

ORDERING/TECHNICAL INFORMATION: PHONE 608-274-4330

www.promega.com • ©2010 PROMEGA CORPORATION • ALL RIGHTS RESERVED • PRICES AND SPECIFICATIONS SUBJECT TO CHANGE WITHOUT PRIOR NOTICE • PRINTED IN USA, REV 5/10 • 18744 • PART #FB087

