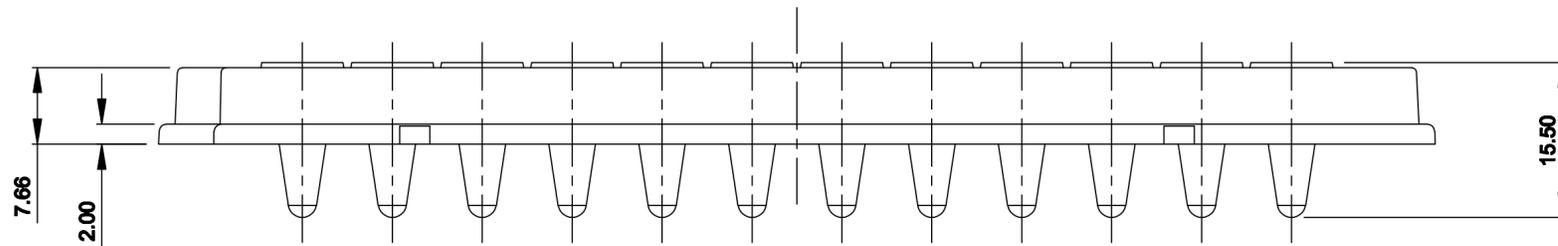
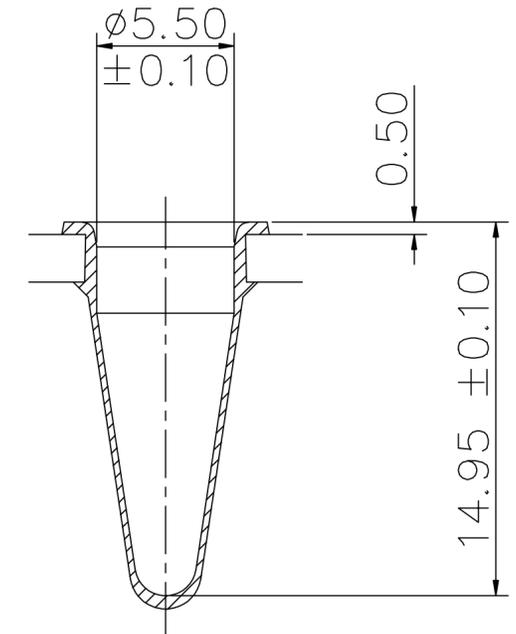
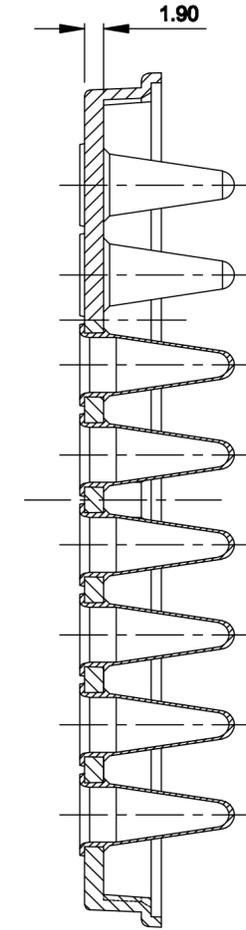
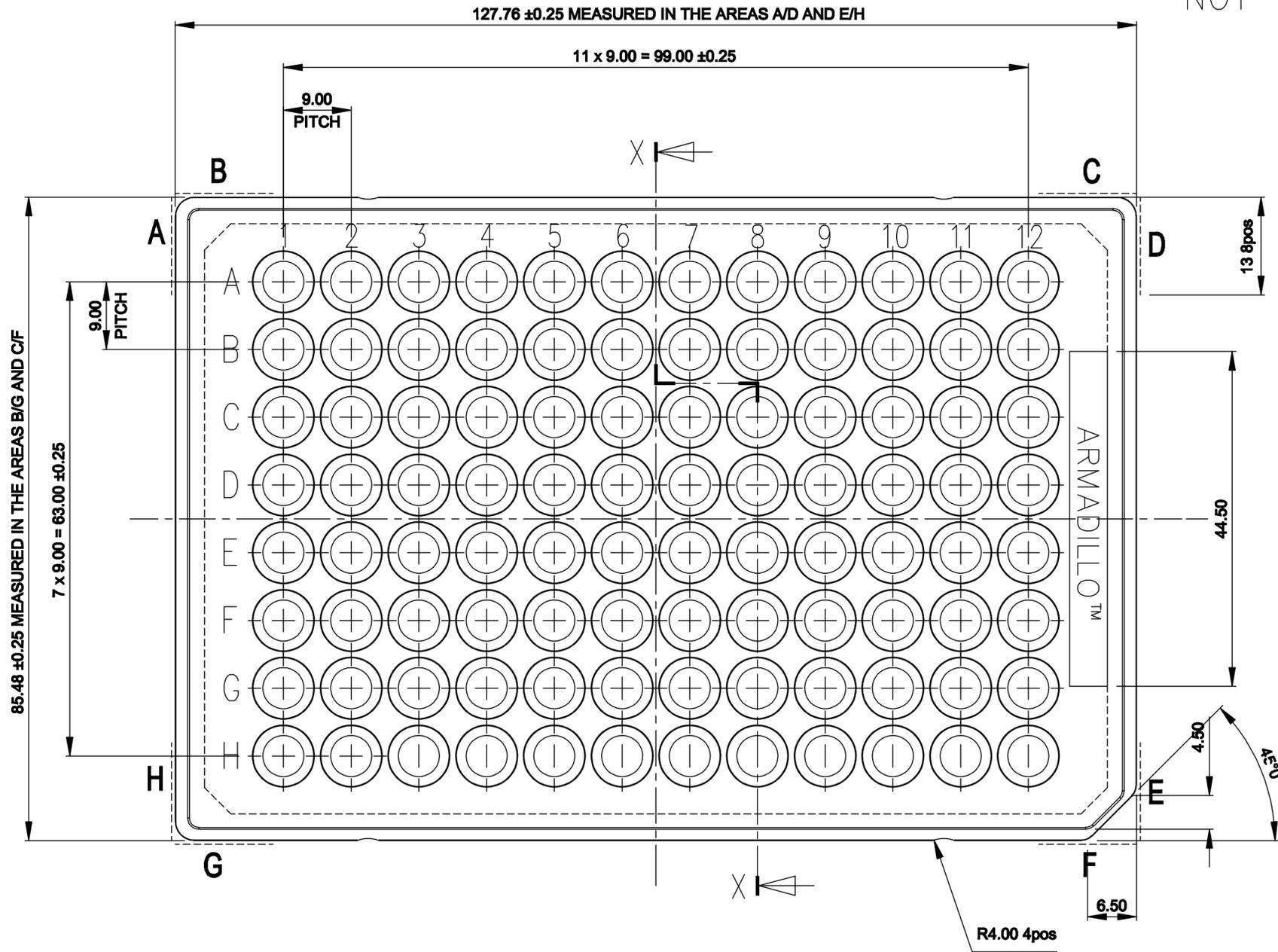


NOT TO SCALE

Copyright Thermo Fisher Scientific  
This drawing and/or any technical information relating to it is issued in strict confidence and may not be reproduced or used for manufacture without written permission. Upon receipt of which it is the users responsibility to seek confirmation of any dimensions and production tolerances.



Pub No. MAN0014463

All dimensions in mm

TITLE		
Semi Skirted Low Profile 96 Well Armadillo PCR Plate		
DRG. NO.	ISSUE	DATE
AB-2496	ORIGINAL	5/8/14