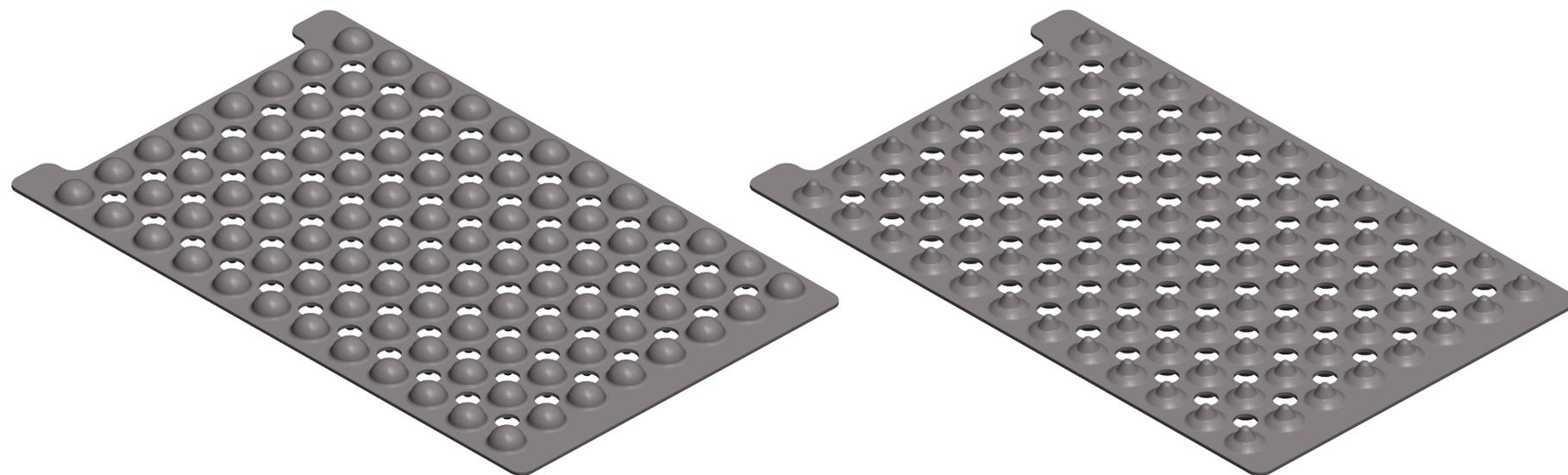


SECTION A-A



PART #	PACK SIZE	DESCRIPTION
N8010550	5 COVERS	MicroAmp® 96-Well Full Plate Cover

<small>UNLESS OTHERWISE SPECIFIED, DIMENSIONS AND TOLERANCES ARE PER ASME Y14.5M AND ARE IN MILLIMETERS. TOLERANCES ARE AS FOLLOWS:</small>	MATERIAL	N/A									
	FINISH	N/A									
	DESIGNED	DATE									
<input checked="" type="checkbox"/>	THIRD ANGLE PROJECTION		<table border="1"> <tr> <td>SIZE</td> <td>SCALE</td> <td>PART NO.</td> <td>REV</td> </tr> <tr> <td>D</td> <td>3.000</td> <td>N8013903</td> <td></td> </tr> </table>	SIZE	SCALE	PART NO.	REV	D	3.000	N8013903	
SIZE	SCALE	PART NO.	REV								
D	3.000	N8013903									

© 2016 ThermoFisher Scientific corporation. All rights reserved. These drawings and specifications are the property of ThermoFisher Scientific and are issued in strict confidence and shall not be reproduced, copied, or used as the basis for manufacture, or sale of, or any unauthorized purposes related to, the matter depicted herein, without the prior written consent of ThermoFisher Scientific corporation.

METRIC

ThermoFisher
SCIENTIFIC

DESCRIPTION
BLANKET-TC11
96 WELL TRAY

DO NOT SCALE DRAWING SHEET 1 OF 1