



PART #	PACK SIZE	DESCRIPTION
4306311	100 FILMS	MicroAmp® Clear Adhesive Film

<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>2.000</td> <td>100065675</td> <td></td> </tr> <tr> <td colspan="2">DO NOT SCALE DRAWING</td> <td colspan="2">SHEET 1 OF 1</td> </tr> </table>	SIZE	SCALE	PART NO.	REV	D	2.000	100065675		DO NOT SCALE DRAWING		SHEET 1 OF 1	
SIZE	SCALE	PART NO.	REV												
D	2.000	100065675													
DO NOT SCALE DRAWING		SHEET 1 OF 1													
<b>METRIC</b>															

© 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.