



SECTION A-A

PART #	PACK SIZE	DESCRIPTION
4375816	20 PLATES	MicroAmp® Fast Optical 48-Well Reaction Plate

<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	
METRIC	THIRD ANGLE PROJECTION		<small>SIZE SCALE PART NO. REV</small> D 3.000 4368897
	<small>DO NOT SCALE DRAWING SHEET 1 OF 1</small>		

© 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.