

D

C

B

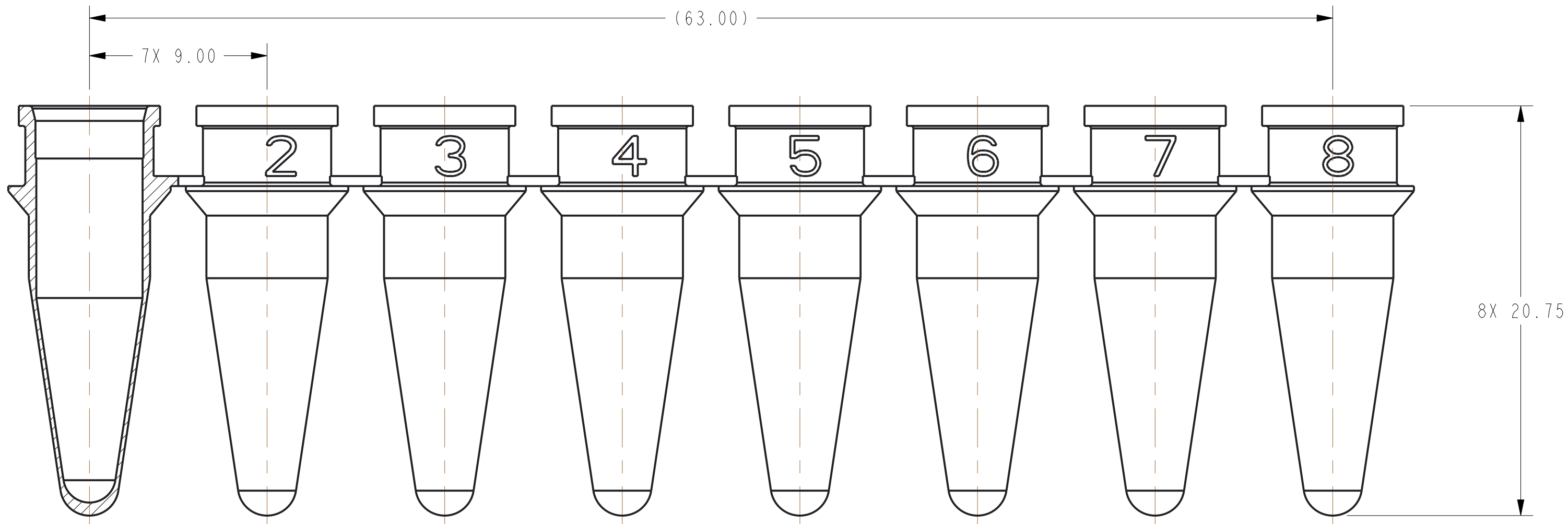
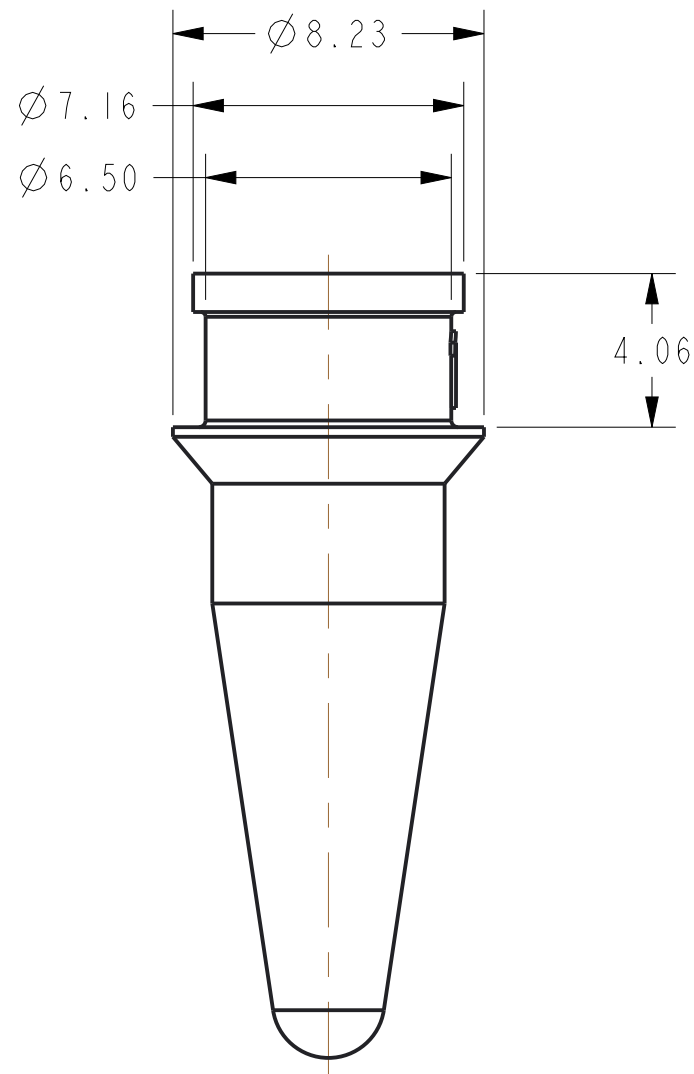
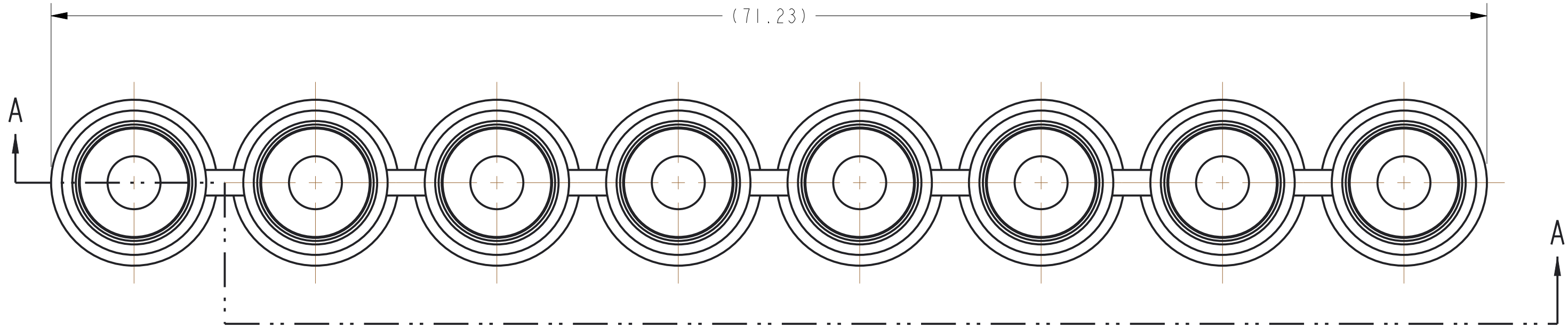
A

D

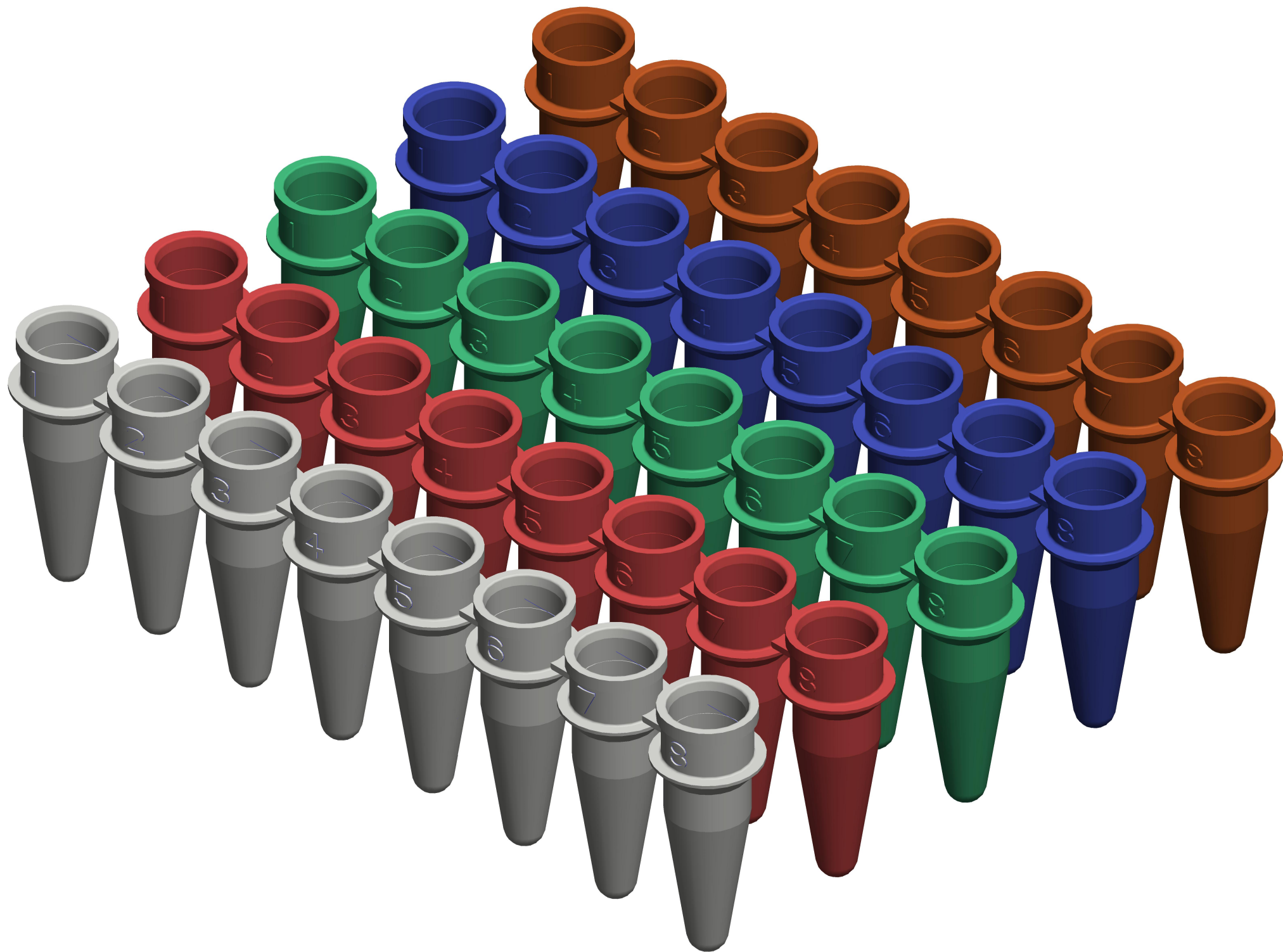
C

B

A




SECTION A-A



PART #	PACK SIZE	DESCRIPTION
N8010580	125 STRIPS	MicroAmp® Optical 8-Tube Strip, 0.2 mL
N8010838	120 STRIPS	MicroAmp® Optical 8-Tube Strip, 0.2 mL, Assorted Colours

© 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 master depicted herein, without the prior written consent of ThermoFisher Scientific corporation.

UNLESS OTHERWISE SPECIFIED, DIMENSIONS AND TOLERANCES ARE PER ASME Y14.5M AND ARE IN MILLIMETERS. TOLERANCES ARE AS FOLLOWS: ✓	MATERIAL N/A		ThermoFisher SCIENTIFIC					
	FINISH N/A							
	DESIGNED		DATE		DESCRIPTION TUBE - TC11 8 / STRING			
	THIRD ANGLE PROJECTION 							
METRIC	SIZE D	SCALE 5.000	PART NO. N8013994	REV				
	DO NOT SCALE DRAWING			SHEET 1 OF 1				