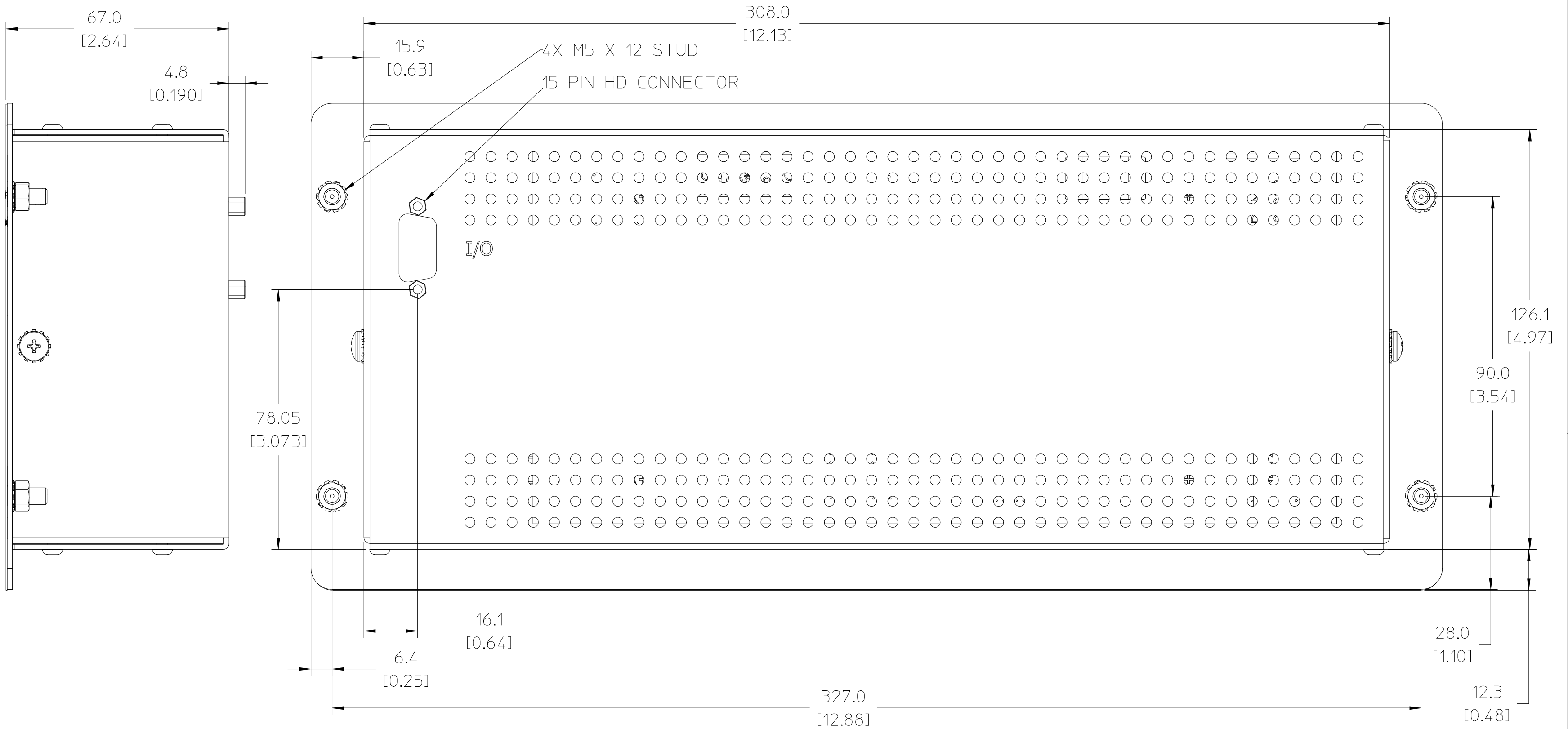
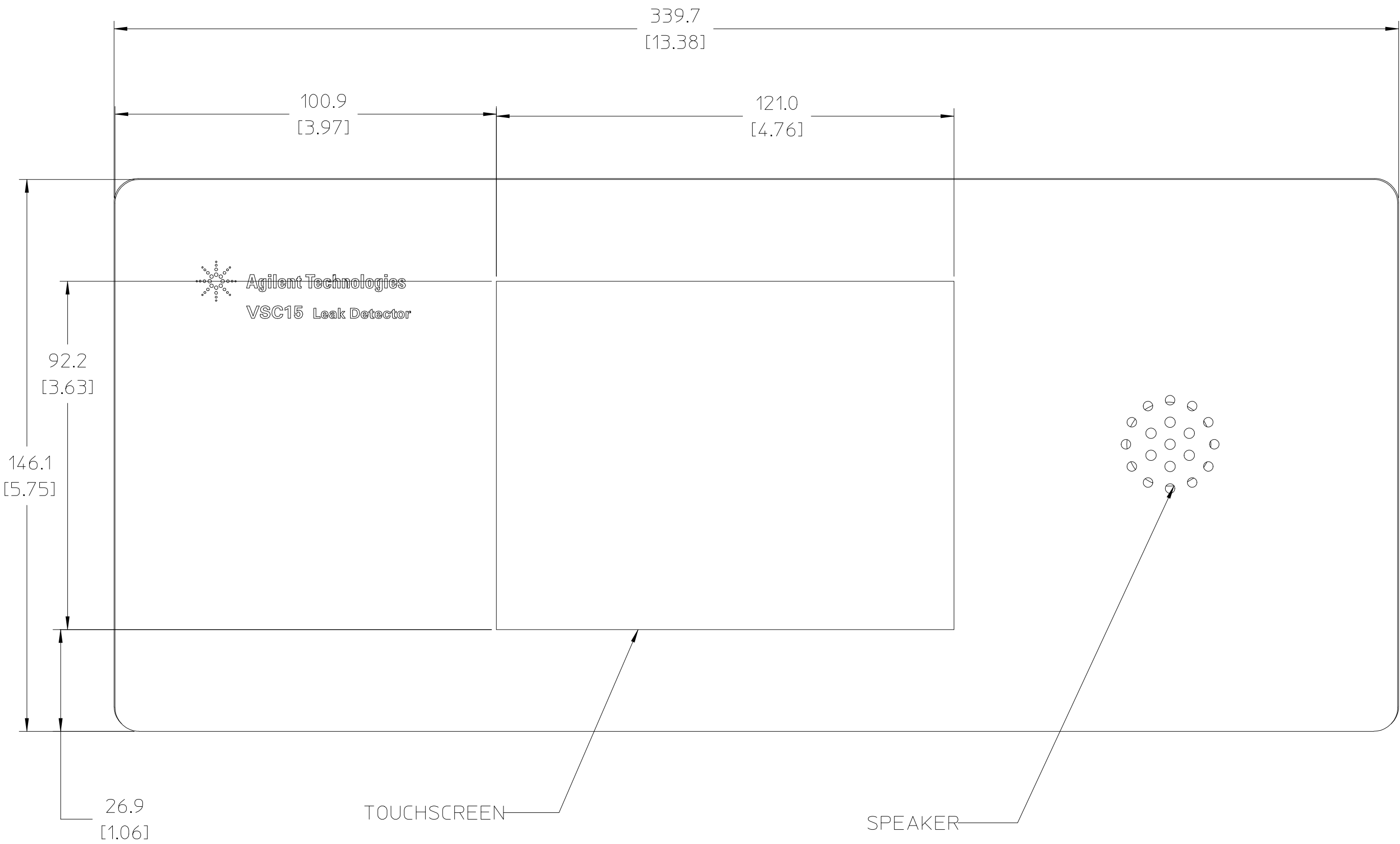




3. ADDITIONAL REQUIREMENTS (PACKAGING, LABELING, QUANTITY, PROCESS REQUIREMENTS, ETC.) ARE SPECIFIED ON PART ORDER.
2. INTERNAL AND EXTERNAL RADII 0.25 mm MAX.
1. FABRICATION TOLERANCES PER A-5951-1561-1 (608) APPLY.

NOTES: (UNLESS OTHERWISE SPECIFIED)