



SECTION DD

Tolerances unless otherwise specified:
 Without Decimal: ± 0.5 mm
 1 Place Decimal: ± 0.1 mm
 2 Place Decimal: ± 0.05 mm
 Material Thickness: ± 0.02 mm
 Undefined Radius: R0.2 mm
 Undefined Draft: 2°
 calculated volume .529 ml

UNLESS OTHERWISE SPECIFIED DIMENSIONS ARE IN MM TOLERANCES ARE: ANGLES: \pm DECIMALS: \pm FRACTIONS: \pm		DWG TITLE		P-98-450V	
APPROVALS	DATE				
DRAWN: CRN	08/22/01				
CHECKED:					
DATE:		SIZE	DWG NO.	REV	
n:\dirge\app\dirge\sl008BAA		D	SL0088	AA	
		SCALE	1 : 1	SHEET 1 OF 1	