



SECTION DD

Tolerances unless otherwise specified:  
 Without Decimal:  $\pm 0.5$  mm  
 1 Place Decimal:  $\pm 0.1$  mm  
 2 Place Decimal:  $\pm 0.05$  mm  
 Material Thickness:  $\pm 0.02$  mm  
 Undefined Radius: R0.2 mm  
 Undefined Draft: 2°  
 Volume of wall: 548 ul

UNLESS OTHERWISE SPECIFIED DIMENSIONS ARE IN MM TOLERANCES ARE: ANGLES: $\pm$ DECIMALS: $\pm$ FRACTIONS: $\pm$		DWG TITLE P-96-450R	
APPROVAL	DATE	SIZE	DWG NO.
DRAWN: CRN	08/22/01	D	SLD087
CHECKED: D.TAUNK		BY	AA
n:\dirga\app\dirg\sl0087AA		SCALE	1 : 1
		SHEET 1 OF 1	