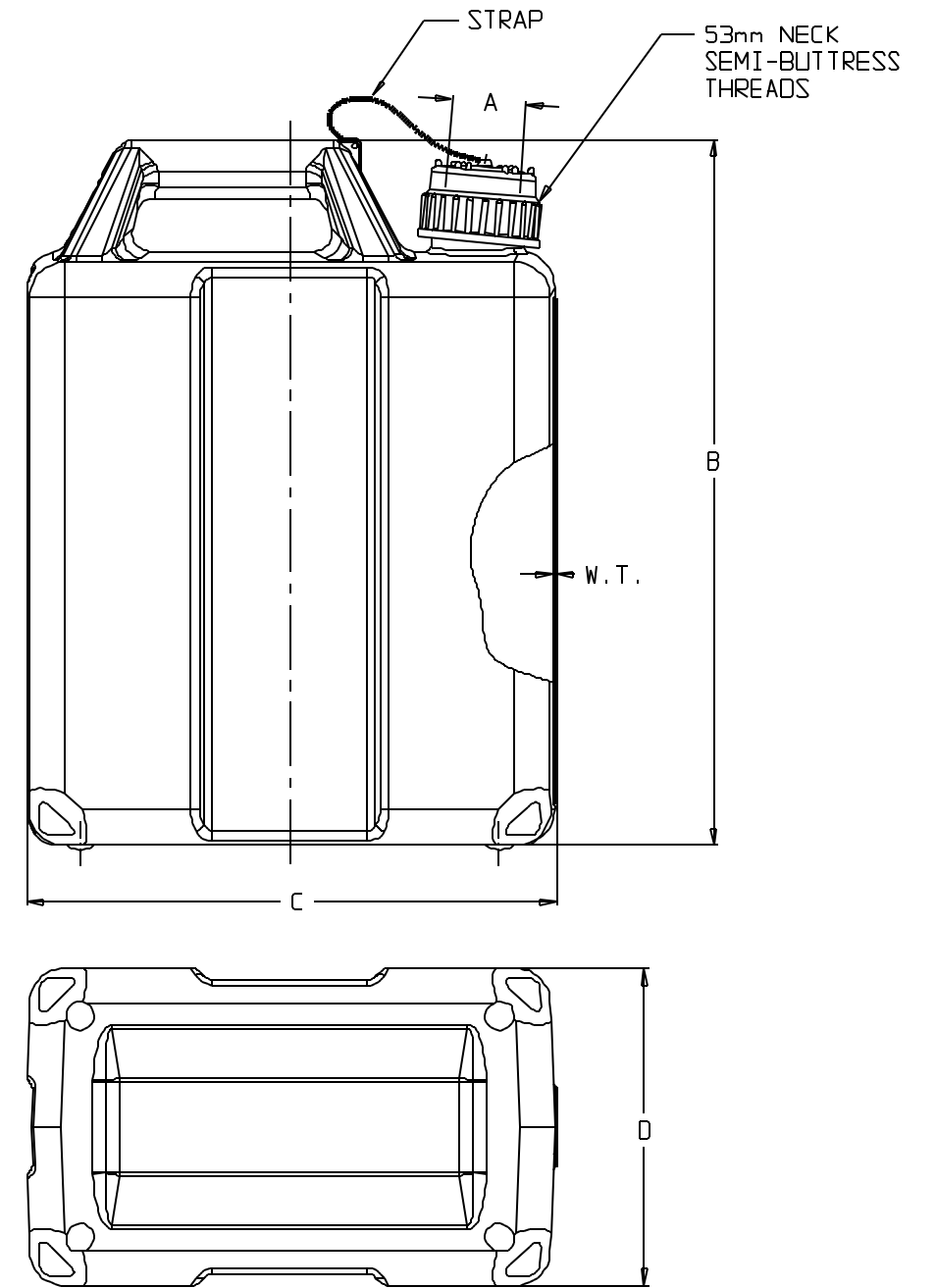


| CATALOG NO. | NOMINAL CAPACITY (L) | MATERIALS | | |
|-------------|-------------------------|-----------|-------------------|-------|
| | | CARBOY | CLOSURE | STRAP |
| 002243-0013 | 13 | HOPE | WHITE PP/NAT HOPE | |

| | | |
|-------------------------------------|--------|-----|
| SIZE (NOM VOL L)..... | 13 | |
| NOM BRIM CAP (L)..... | 14.2 | |
| NECK FINISH (mm)..... | 53 | |
| THREAD (T.P.I.)..... | 4 | |
| NOM. WEIGHT W/ CLOSURE(gms)..... | 1250 | |
| | INCHES | MM |
| A (NECK I.D. AT TOP)..: | 1.5 | 38 |
| B (OVERALL HEIGHT)..: | 14.9 | 378 |
| C (WIDTH)..... | 11.4 | 290 |
| D (DEPTH)..... | 7.45 | 189 |
| W.T.(NOM WALL THK)..: | 0.1 | 2.5 |

2. This is not a production drawing, but is for informational purposes only, and is subject to change without notice.

1. All dimensions are nominal and in inches and millimeters unless otherwise noted.



SPECS ON BACK

NOT TO SCALE

CONFIDENTIAL

THIS PRODUCT SPECIFICATION SHEET CONTAINS CONFIDENTIAL AND/OR PROPRIETARY INFORMATION AND IS TO BE USED ONLY FOR THE EXPRESSED PURPOSES AGREED UPON BY AN AUTHORIZED OFFICER OF THERMO FISHER SCIENTIFIC.

ThermoFisher
SCIENTIFIC

Thermo Fisher Scientific
75 Ponorona Creek Drive
Rochester, NY 14625-2385, USA

| | | | | |
|-------------------|-----|-----------|------------------|------|
| MADE | PB | APP'D PVC | PSS 00-2243-0013 | REV. |
| CHK'D | PVC | 12/4/01 | | 1201 |
| DESCRIPTION: | | | | |
| 13-LITER JERRICAN | | | | |