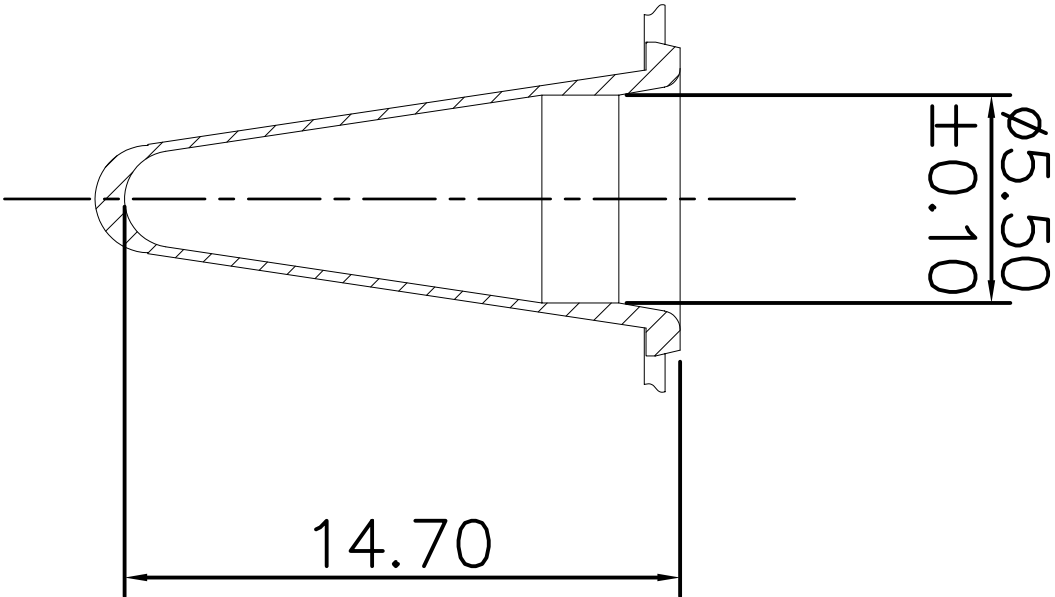
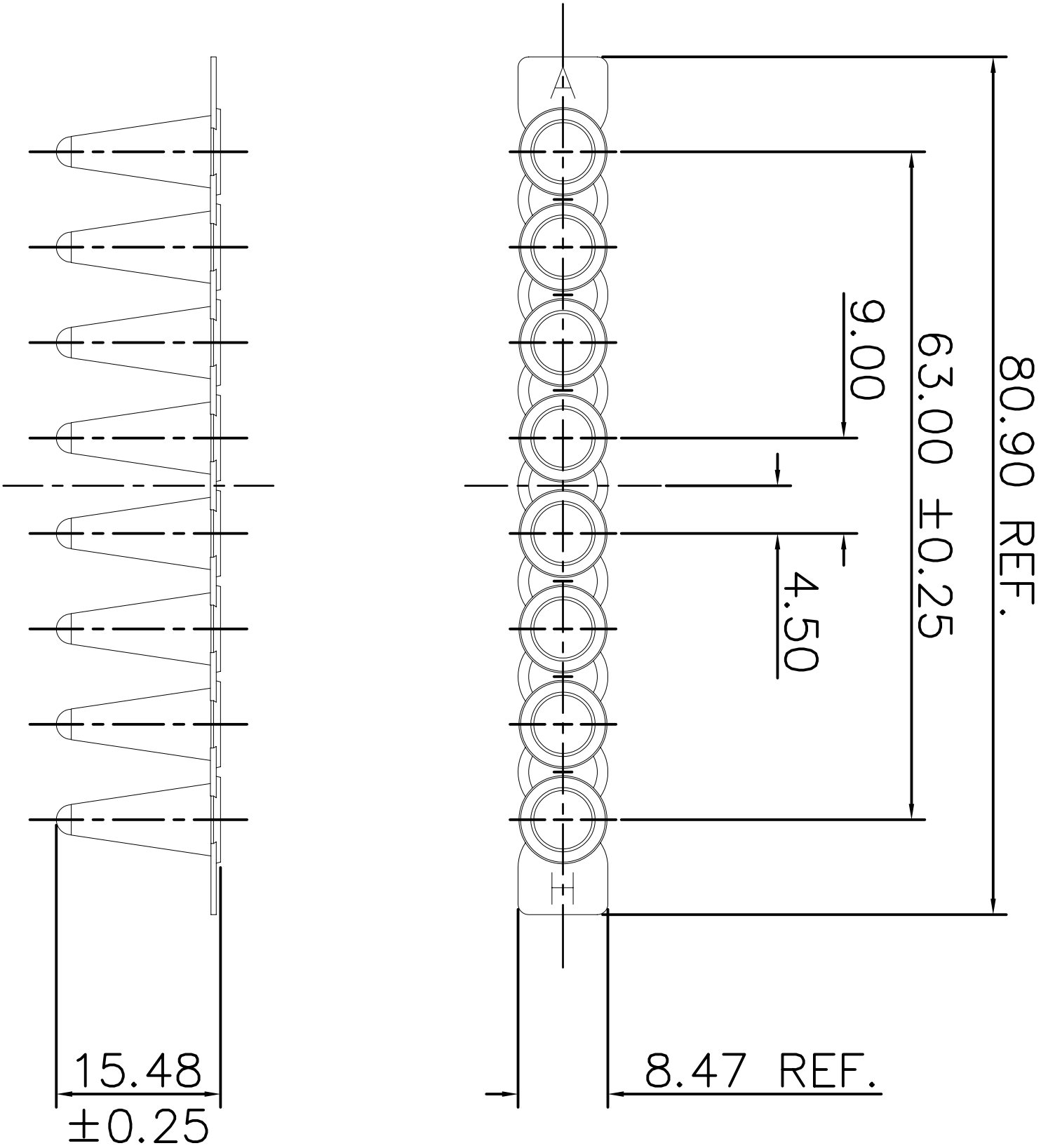


NOT TO SCALE

Copyright ThermoFisher Scientific  
This drawing and/or any technical information relating to it is issued in strict confidence and may not be reproduced or used for manufacture without written permission. Upon receipt of which it is the users responsibility to seek confirmation of any dimensions and production tolerances.



VIAL DETAIL 8pos

All dimensions in mm

| TITLE                        |          |         |  |
|------------------------------|----------|---------|--|
| Strip of 8 Low Profile Tubes |          |         |  |
| DRG. NO.                     | ISSUE    | DATE    |  |
| AB-0771                      | ORIGINAL | 2/11/06 |  |