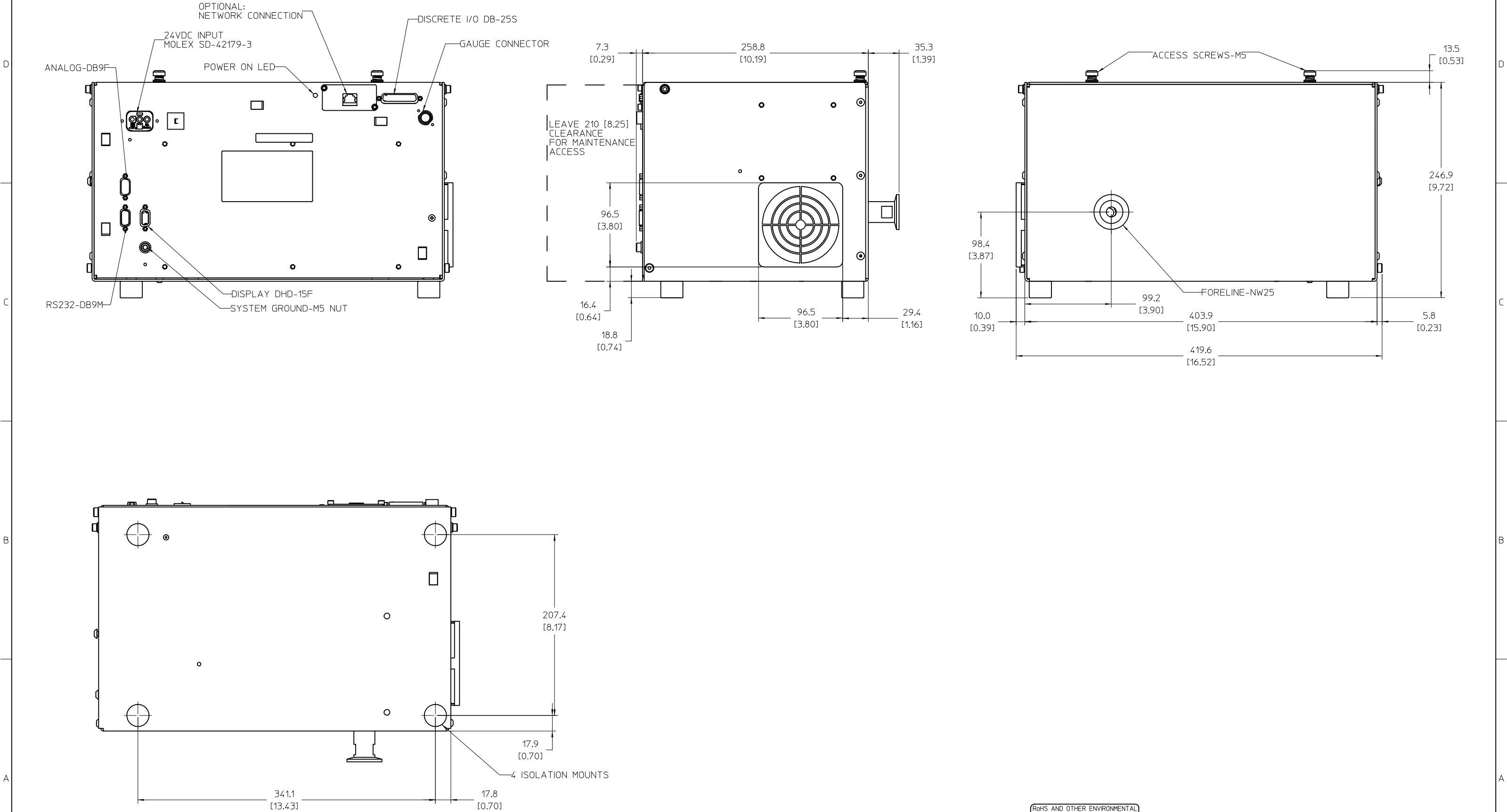


AGILENT CONFIDENTIAL  
THIS DOCUMENT AND ANY ASSOCIATED DATA  
CONTAIN CONFIDENTIAL INFORMATION THAT IS  
AGILENT TECHNOLOGIES PROPERTY. ONLY  
DISCLOSE OR DUPLICATE FOR OTHERS AS  
AUTHORIZED BY AGILENT TECHNOLOGIES.

ASSOCIATED 3D MODEL IS  
GEOMETRY MASTER  
DO NOT MODIFY THE GEOMETRY IN THIS  
2D DRAWING. GEOMETRY MODIFICATIONS  
MUST BE MADE TO THE MASTER 3D MODEL  
AND USED TO UPDATE THIS DRAWING.

| FILE NAME             |  | REVISION | STORE DATE | DOCUMENT NUMBER |            |
|-----------------------|--|----------|------------|-----------------|------------|
| * PER MODEL MANAGER * |  | B        | 2020-01-03 | R3355301        |            |
| SYN                   |  |          | REVISIONS  |                 | REVISED BY |
| B                     |  |          | AS ISSUED  |                 | D. WALL    |
|                       |  |          |            |                 | DATE       |
|                       |  |          |            |                 | 2020-01-03 |



1. WEIGHT 20LBS

NOTES: (UNLESS OTHERWISE SPECIFIED)

|                                                               |  |                                                                                                                |  |                   |  |                                       |  |                                                                                                                                                                       |  |                                |  |
|---------------------------------------------------------------|--|----------------------------------------------------------------------------------------------------------------|--|-------------------|--|---------------------------------------|--|-----------------------------------------------------------------------------------------------------------------------------------------------------------------------|--|--------------------------------|--|
| RoHS AND OTHER ENVIRONMENTAL<br>REQUIREMENTS PER 5951-1745-2. |  |                                                                                                                |  |                   |  |                                       |  |                                                                                                                                                                       |  |                                |  |
| STANDARD TOLERANCES<br>(UNLESS OTHERWISE SPECIFIED)           |  | D. WALL                                                                                                        |  | 2020-01-03        |  | INTERFACE DRAWING,<br>VSC15 BASE UNIT |  |  <b>Agilent Technologies</b><br><small>© Agilent Technologies, Inc. 2016</small> |  | PART<br>NUMBER<br><br>R3215301 |  |
| LINEAR:                                                       |  | DRAWN BY                                                                                                       |  | DATE (YYYY-MM-DD) |  |                                       |  |                                                                                                                                                                       |  |                                |  |
| ANGULAR:                                                      |  | NAME                                                                                                           |  | 2020-01-03        |  |                                       |  |                                                                                                                                                                       |  |                                |  |
|                                                               |  | ENGINEER/CHECKER                                                                                               |  | DATE              |  |                                       |  |                                                                                                                                                                       |  |                                |  |
|                                                               |  | UNLESS OTHERWISE SPECIFIED                                                                                     |  |                   |  |                                       |  |                                                                                                                                                                       |  |                                |  |
|                                                               |  | * DIMENSIONS ARE IN MILLIMETERS                                                                                |  |                   |  |                                       |  |                                                                                                                                                                       |  |                                |  |
|                                                               |  | * THIRD ANGLE PROJECTION  |  |                   |  |                                       |  |                                                                                                                                                                       |  |                                |  |
|                                                               |  | * ADHERENCE TO DWG A-5951-6561-1<br>(COMPANY STD. SECTION 608) AND<br>A-5951-1745-1 (GSE) IS REQUIRED          |  |                   |  |                                       |  |                                                                                                                                                                       |  |                                |  |
| DO NOT SCALE THIS                                             |  | GD&T SYMBOLOGY<br>PER                                                                                          |  | TITLE             |  | FRAME DATE: 2021-01-01                |  |                                                                                                                                                                       |  |                                |  |
| METRIC<br>DRAWING                                             |  | ASME Y14.5-2009                                                                                                |  | 1 OF 1            |  | 0.5X                                  |  | D                                                                                                                                                                     |  | R3355301                       |  |
|                                                               |  |                                                                                                                |  | SHEET             |  | SCALE                                 |  | SIZE                                                                                                                                                                  |  | DOCUMENT NUMBER                |  |