

## TDS OptiPlate-24 / CulturPlate-24 (PN 6005186 / 6005168)

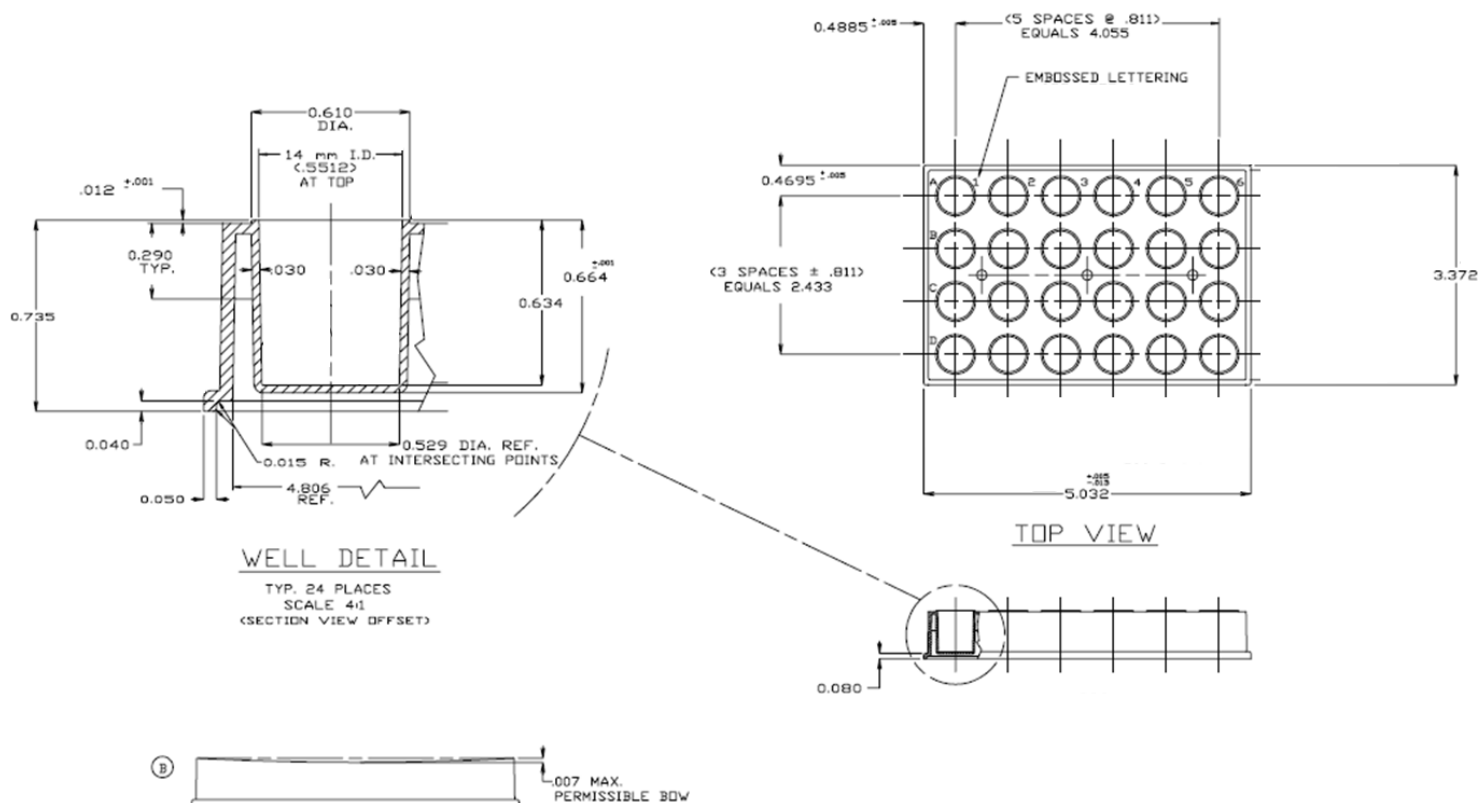
### Material and surface treatment

Opaque white 24-well microplate made of polystyrene. OptiPlate-24 (6005186) is made of untreated polystyrene and CulturPlate-24 (600516) is sterilized and tissue culture treated to facilitate cell adherence and proliferation.

### Packaging

6005186 is a box with 100 plates packed as 4 bags with 25 plates each. 6005168 is a case with 50 individually pouched, lidded microplates.

### Dimensions --- customer drawing



All dimensions in inches

Product distributed by:  
PerkinElmer Health Sciences B.V.  
The Netherlands