

Research use only. Not for use in diagnostic procedures.

Technical datasheet:

Dimensions of Low-Profile Microplate Lid for PhenoPlate Imaging Microplates

**Part numbers:** PN 6000031 Clear, sterile, low profile, for 'PhenoPlate' imaging

**Material and treatment:** Clear, low profile, polystyrene microplate lid. Fits on all imaging microplates of the 'PhenoPlate' - range.

**Packaging:** 6000031 is a case with 200 lids.

Dimensions -- Customer Drawing:

