

Research use only. Not for use in diagnostic procedures.

Technical datasheet: Dimensions of Clear Universal Microplate Lid

Part numbers: PN 6000030 Clear, untreated

Material and treatment: Clear, automation-friendly polystyrene microplate lid. Fits on all microplates and (polypropylene) reservoirs. Lid is available in 'untreated' version only.

Packaging: 6000030 is a case with 100 lids, packed as 5 bags of 20 lids each.

Dimensions -- Customer Drawing:

