

PRODUCT NAME

96 well microplates

Cat. No: **WA-003-0010**

Format: 10 pc

Description:

The 96 well microplates with U-bottom offered by Diagenode are made by polystyrene and are compatible with automated systems.

The product is offered in clear colour.

Plates are free of heavy metals and are free of detectable DNase/RNase, human DNA and pyrogens.

Working temperatures are from -20° to 60° degrees.

Storage is recommended at room temperature.

Autoclavability is not possible and centrifugation must be done at a maximal of 4500xg in a swinging-bucket rotor

Dimensions:

Length: 127.76 mm (+/- 0.2mm)

Width: 85.48 mm (+/- 0.2 mm)

Curvature: ~200 µm

Volume per well:

Total volume: 320 µl

Working volume: 40 to 280 µl