

Research use only. Not for use in diagnostic procedures.

Ergonomic High-throughput Cell Counting Plates, 12 x 2 Orientation (CHM24-A100)

Material: Clear polystyrene all-plastic plate.

Packaging:

Part Numbers	Format	Number of Plates
CHM24-A100-098	One sleeve of 10 plates	10
CHM24-A100-099	One case containing 4 sleeves, 10 plates per sleeve	40
CHM24-A100-100	Five cases, each case containing 4 sleeves, 10 plates per sleeve	200

Dimensions – Customer Drawing:

Drawing is for reference only. All dimensions are in millimeters (mm).

