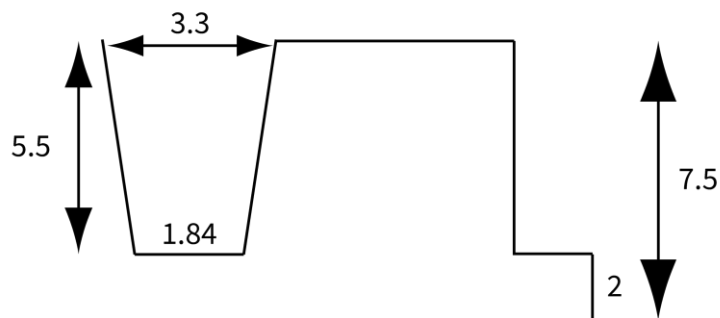
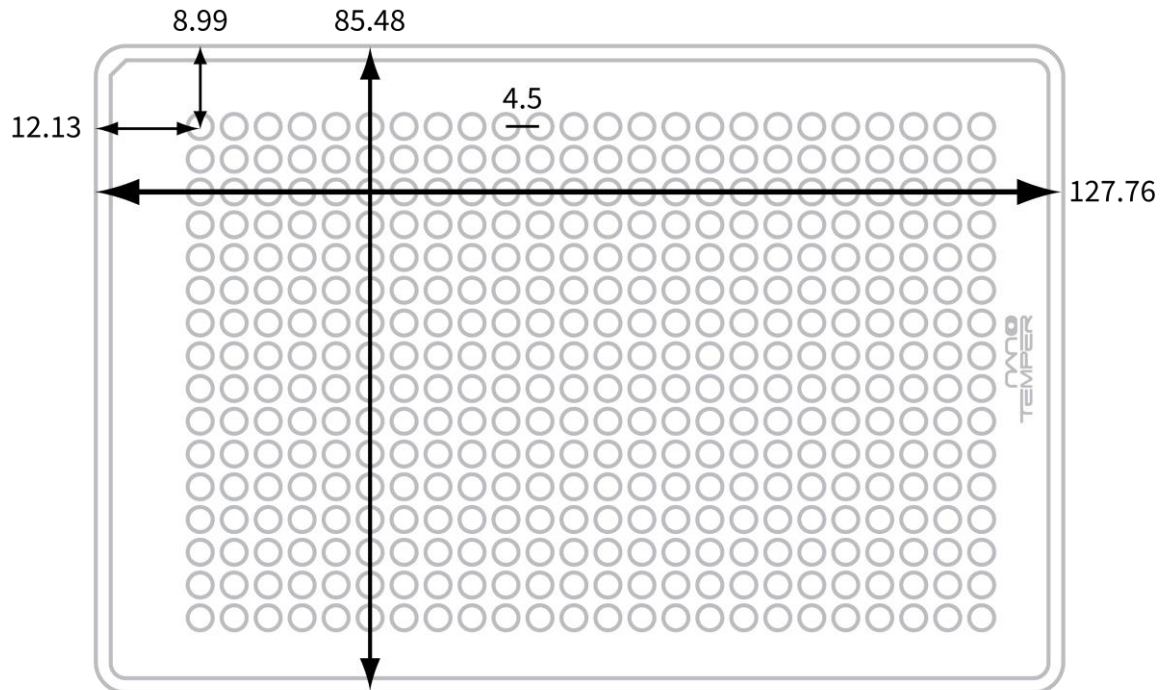


Dianthus Plate Specifications

Measures (in mm):



Important information:

Working volume: 18-25 μ L

Coated surface to avoid protein or nucleic acid adsorption

Conical well geometry

Temperature tolerance: -20°C – +60°C