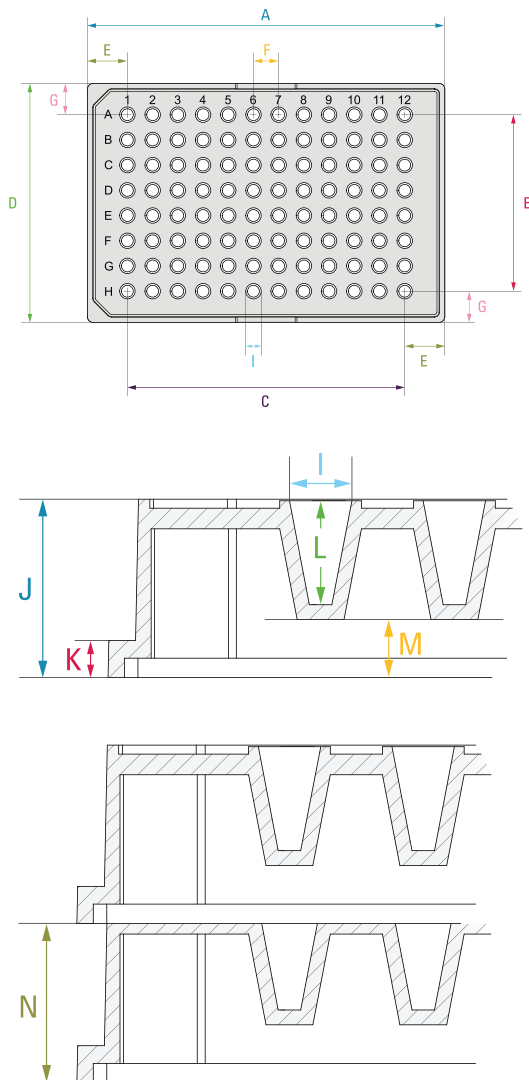


Research use only. Not for use in diagnostic procedures.

# HTRF 96-well low volume

## 96-Well plate

Product	Part Number	Well Density	Color	Treatment	Quantity
HTRF Low Volume Plate	66PL96001	96	White	Untreated	1
	66PL96005	96	White	Untreated	5
	66PL96025	96	White	Untreated	25
	66PL96100	96	White	Untreated	100



Parameter	Specification
Material	Polystyrene
Well Density	96-well
Well Shape	Round
Well-bottom Shape	V-bottom
A. Plate Length (mm)	127.6
B. Row Coordinate (mm)	63.0
C. Column Coordinate (mm)	99.0
D. Plate Width (mm)	85.4
E. A1 Column Offset (mm)	14.3
F. Well to Well Spacing (mm)	9.0
G. A1 Row Offset (mm)	11.2
I. Well Diameter (mm)	4.2
J. Plate Height (mm)	12.0
K. Flange/Skirt/Border height (mm)	2.5
L. Well Depth (mm)	7.0
M. Well Bottom Elevation (mm)	4.0
N. Stack Height (mm)	10.5
Maximum Well Volume	40µL
Working Volume	20µL