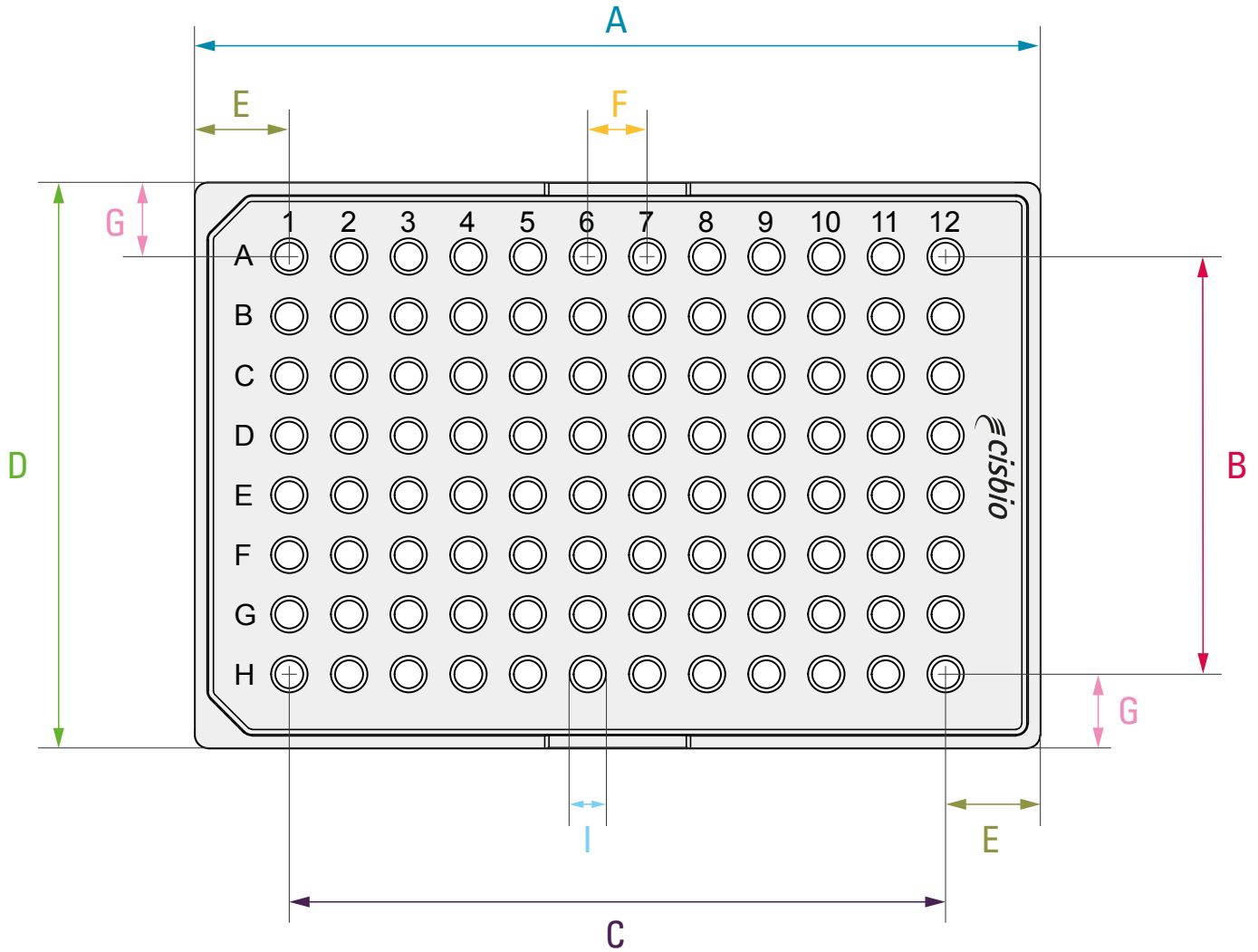


HTRF 96-WELL LOW VOLUME WHITE PLATE CHARACTERISTICS

TECHNICAL INFORMATION



A = 127.60 mm (Plate length)

B = 63 mm (Row coordinate of bottom left or right corner)

C = 99 mm (Column coordinate of top or bottom right corner well)

D = 85.40 mm (Plate width)

E = 14.3 mm (Column offset coordinate of top - or bottom left corner well)

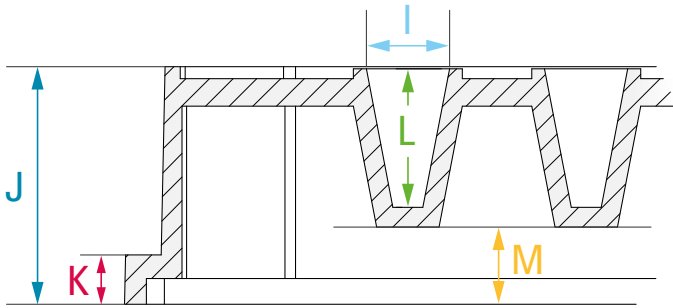
F = 9 mm Well spacing (center to center)

G = 11.2 mm (Row offset coordinate of top or bottom left corner well)

I = 4.20 mm (Well diameter top)

D - G (bottom right Y) = 74.2mm
A - E (bottom right X) = 113.3mm

Working volume = 20µL
Maximum volume = 40µL



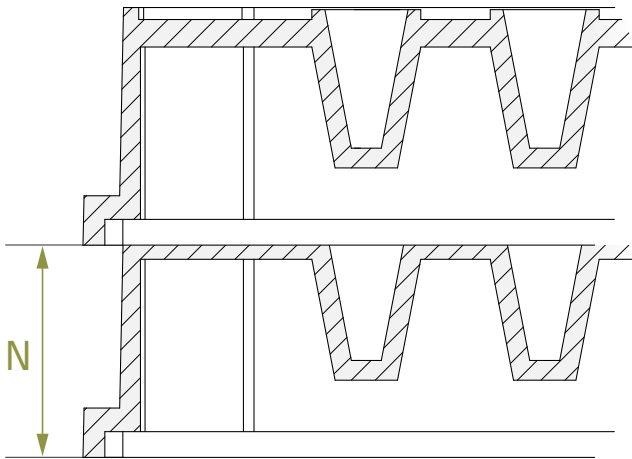
I = 4.20 mm (Well diameter top)

J = 12 mm (Plate height)

K = 2.50 mm (Border)

L = 7 mm (Well depth)

M = 4 mm (Well bottom elevation)



N = 10.50 mm (Stack height)

HTRF® 96-WELL PLATE

Product	Size	Part #
HTRF 96 well low volume plate	1 plate	66PL96001
	5 plates	66PL96005
	25 plates	66PL96025
	100 plates	66PL96100

For more information, visit
www.cisbio.com/HTRF-96-well-plate



FOR MORE INFORMATION

Europe and other countries +33(0)466 796 705 U.S. and Canada 1-888-963-4567 China +86 10 8080 9288

Visit www.htrf.com to find a list of our regional distributors