
Streptavidin-Coated Strip Microwell Plate (Flat)

(Catalog No. R-1102)

Description

Streptavidin-coated surfaces offer a powerful and universal method to bind any biotinylated molecule (proteins, peptides, polysaccharides, oligonucleotides, DNA fragments, etc.). Epigentek's Streptavidin-Coated Strip Microwell Plate has the following features:

- Perforated 8-well strips which fit onto both 8x12 and single well holding frames
- Pre-blocked and ready-to-use
- High sensitivity to detect as low as 0.05 ng/well of biotinylated-HRP with a detection range of 0.05 to 10 ng/well of biotinylated molecules
- SBS standard compliant dimensions
- Ideal for fast processing on automated machinery and instruments

Storage Temp & Conditions

Store aliquots at 4°C.

Application Notes

Used for the detection of various concentration of biotinylated molecules and general ELISA screening

Ordering Information:

Products	Size	Cat. No.
Streptavidin-Coated Strip Microwell Plate (Flat)	5/pk	R-1102 -1

This product is for research purposes only. Not intended for use in diagnostic procedures.