
Streptavidin

(Catalog No. R-1100)

Description

Streptavidin is a tetrameric protein, with each subunit able to bind a single molecule of the vitamin biotin. Streptavidin-biotin bond is the strongest known non-covalent interaction with $K_d \sim 10^{-15}$ M, and exhibits less nonspecific binding than avidin. Streptavidin is used extensively in molecular biology and bionanotechnology due to the streptavidin-biotin complex's resistance to organic solvents, denaturants, detergents, proteolytic enzymes, and extremes of temperature and pH.

Molecular Weight (kDa)

60 kDa

Formulation

10 mM phosphate buffered saline, pH 7.4

Source

Expressed in E coli

Size

1 mg

Storage Temp & Conditions

>6 months at -20°C . Avoid multiple freeze/thaw cycles.

Application Notes

Useful for the purification or detection of various biomolecules

Ordering Information:

Products	Size	Cat. No.
Streptavidin	1 mg	R-1100 -1

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