

Streptavidin, Recombinant

Catalog Number: AS-72177-100

Amount: 100 mg

Physical State: Lyophilized powder.

Storage: The lyophilized powder is stable for years when stored at 4°C with

desiccant. Reconstitute the powder with desired buffer. The

reconstituted solution will be stable for at least six months at -20°C The reconstituted solution may also be divided into aliquots to

avoid repeated freeze-thaw cycles.

Instruction: Streptavidin is a nonglycosylated tetrameric protein that binds

biotin noncovalently and with high affinity (Kd $\sim 10^{-15}$ M). The recombinant streptavidin was expressed in *E. coli*. It shows one

major band about 56 kDa on SDS-PAGE. The activity of recombinant streptavidin is ~17 units/mg, where one unit is defined as the amount of protein required to bind 1.0 μ g of D-biotin. The extinction coefficient of streptavidin is $E^{1\%}$ 280=32;

1 mg/mL $A_{280} = 3.2$.

For *in vitro* research use only This product not of animal origin