



Streptavidin, Recombinant

Catalog Number: AS-72177-5

Amount: 5 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 ($K_d \sim 10^{-15}$ 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 µg of D-biotin. The extinction coefficient of streptavidin is $E_{280}^{1\%} = 32$; 1 mg/mL $A_{280} = 3.2$.

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