

## Streptavidin, Recombinant

Catalog Number: AS-72177-500

**Amount:** 500 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  $\mu$ 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