

Streptavidin, Recombinant

Catalog Number:	AS-72177-1000
Amount:	1 g
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 ~10 ⁻¹⁵ M). The recombinant streptavidin was expressed in <i>E. coli</i> . 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