

Product Information Sheet

Product Name: Streptavidin

Catalog Number: AS-60659

Amount: 5 mg (This product containing 90% streptavidin and 10% sodium

chloride by weight.)

Physical State: Lyophilized powder.

Storage: The lyophilized powder is stable for years at 4 °C. Reconstitute the

powder with desired buffer, such as phosphate-buffered saline. The reconstituted solution will be stable for three months at 4°C if it is added with sodium azide to a final concentration of 2 mM or thimerosal to 0.02%. The reconstituted solution may also be divided into aliquots and stored at -20°C for long-term storage.

Description: Streptavidin is a 55,000-dalton nonglycosylated tetrameric protein.

It was purified from culture broth of the bacterium *Streptomyces avidinii*. Streptavidin binds to D-biotin noncovalently and with high affinity ($K_d \sim 10^{-15} M$). The activity of this product is ~ 14 units/mg, where one unit is defined as the amount of protein required to bind 1.0 µg of D-biotin. The concentration (mg/mL) of

streptavidin is calculated as A_{280nm} / 3.2.

For in vitro research use only