
**Blotting Grade
Streptavidin
(from *S. avidinii*)**

**Catalog Number
170-6408**



Specifications

Contents 1 mg

Storage Store at -20 °C. As a lyophilized powder, streptavidin is stable for 1 year at -20 °C. Upon rehydration in buffer or deionized water, aliquot and store at -20 °C. Avoid repeated freeze/thaw cycles. Rehydrated streptavidin is stable for at least 6 months at -20 °C.

Note: For research use only. Not for drug or human use.