

Biotin & Streptavidin Products

Limited Time Only

**25%
OFF**

Use promo code **SB25**

Coupon code must be provided at the time of sale for promotion to apply. Offer only applies to purchases of biotin & streptavidin products. Unless otherwise stated, offer only applies to purchases directly from AAT Bioquest, Inc. Contact us for purchases outside of USA. Offer expires on 07/31/2019.