



Product Information Sheet

Product Name: Streptavidin, Horseradish Peroxidase Conjugated

Catalog Number: 60668

Amount: 1 mg

Concentration: 1 mg/mL

Degree of Substitution: HRP/protein mole ratio is labeled on each vial.
(DOS)

Buffer: 10 mM phosphate, 150 mM NaCl, pH 7.2, with 0.5% Proclin 300 as preservative.

Storage: Store at 4°C. DO NOT FREEZE.

Instructions: The suggested concentration for Western blot, ELISA, and immunohistochemistry is 0.1-1 µg/mL.

For *in vitro* research use only