



43-4301 1 mg
Streptavidin
(*S.avidinii*)
Lot No.

STREPTAVIDIN (from *S.avidinii*)

FORM:

Streptavidin is provided in lyophilized form in 10 mM phosphate buffered saline, pH 7.4. Reconstitute with 1 ml of distilled or deionized water.

CAPACITY:

Binding activity of streptavidin is determined to be ≥ 11 units per mg protein. One unit of binding activity is defined as that amount of protein which binds 1 μ g of biotin (Green, N.M., Biochem. J., 89:599,1963)

STORAGE:

Store at 4°C.

REMARKS:

For research use only. Not for human or drug use. No preservative added.

www.invitrogen.com

Invitrogen Corporation • 542 Flynn Rd • Camarillo • CA 93012 • Tel: 800.955.6288

E-mail: techsupport@invitrogen.com

PI 43-4301

(Rev 09/08) DCC-08-1572

Important Licensing Information - These products may be covered by one or more Limited Use Label Licenses (see the Invitrogen Catalog or our website, www.invitrogen.com). By use of these products you accept the terms and conditions of all applicable Limited Use Label Licenses. Unless otherwise indicated, these products are for research use only and are not intended for human or animal diagnostic, therapeutic or commercial use.



43-4301 1 mg
Streptavidin
(*S.avidinii*)
Lot No.

STREPTAVIDIN (from *S.avidinii*)

FORM:

Streptavidin is provided in lyophilized form in 10 mM phosphate buffered saline, pH 7.4. Reconstitute with 1 ml of distilled or deionized water.

CAPACITY:

Binding activity of streptavidin is determined to be ≥ 11 units per mg protein. One unit of binding activity is defined as that amount of protein which binds 1 μ g of biotin (Green, N.M., Biochem. J., 89:599,1963)

STORAGE:

Store at 4°C.

REMARKS:

For research use only. Not for human or drug use. No preservative added.

www.invitrogen.com

Invitrogen Corporation • 542 Flynn Rd • Camarillo • CA 93012 • Tel: 800.955.6288

E-mail: techsupport@invitrogen.com

PI 43-4301

(Rev 09/08) DCC-08-1572

Important Licensing Information - These products may be covered by one or more Limited Use Label Licenses (see the Invitrogen Catalog or our website, www.invitrogen.com). By use of these products you accept the terms and conditions of all applicable Limited Use Label Licenses. Unless otherwise indicated, these products are for research use only and are not intended for human or animal diagnostic, therapeutic or commercial use.