



43-4314 1 mg
TRITC-STREPTAVIDIN

TRITC-STREPTAVIDIN

FORM:

1 ml liquid. TRITC-Streptavidin is provided in 10 mM phosphate-buffered saline, pH 7.4, containing 1% bovine serum albumin, 40% glycerol, and 0.1% sodium azide.

Streptavidin (from *S. Avidinii*) is conjugated with TRITC under optimum condition. The conjugate is purified by molecular sieving and is completely free of unbound TRITC.

APPLICATIONS:

TRITC-Streptavidin can be used for the detection or localization of biotinylated biomolecules. Working dilution for specific application should be determined by the investigator to obtain the best conditions.

STORAGE:

Store at 4°C. Avoid exposure to light.

REMARKS:

For research use only. Not for human or drug use. Product contains 0.1% sodium azide as preservative.

www.invitrogen.com

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

E-mail: techsupport@invitrogen.com

PI434314

(Rev 09/09) DCC-08-1089

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.



43-4314 1 mg
TRITC-STREPTAVIDIN

TRITC-STREPTAVIDIN

FORM:

1 ml liquid. TRITC-Streptavidin is provided in 10 mM phosphate-buffered saline, pH 7.4, containing 1% bovine serum albumin, 40% glycerol, and 0.1% sodium azide.

Streptavidin (from *S. Avidinii*) is conjugated with TRITC under optimum condition. The conjugate is purified by molecular sieving and is completely free of unbound TRITC.

APPLICATIONS:

TRITC-Streptavidin can be used for the detection or localization of biotinylated biomolecules. Working dilution for specific application should be determined by the investigator to obtain the best conditions.

STORAGE:

Store at 4°C. Avoid exposure to light.

REMARKS:

For research use only. Not for human or drug use. Product contains 0.1% sodium azide as preservative.

www.invitrogen.com

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

E-mail: techsupport@invitrogen.com

PI434314

(Rev 09/09) DCC-08-1089

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.