

TREVIGEN[®] Product Data

For Research Use Only. Not For Use In Diagnostic Procedures.

Strep-AMCA

Catalog #: 4887-100-03

Size: 100 µl at 1 mg/ml

Description: AMCA-conjugated Streptavidin

Fluorophore: 7-Amino-4-methylcoumarin-3-acetic acid

$A_{\text{max}} = 350 \text{ nm}$

$E_{\text{max}} = 450 \text{ nm}$

Physical State: This conjugate is provided in 50% glycerol.

Preservative: 0.01% Thimerosal, 0.05% Sodium Azide

Storage: Store at -20°C.

Country of Origin: USA

Applications:

Use between 2-5 µg/ml for most applications.

For flow cytometry applications, the conjugate has been used at 2 µg/ml dilution.

Optimization may be required.

Lot Specific Data:

Lot #:

Fluorophore/Protein ratio: $A_{350}/A_{280} = 0.6$

Streptavidin Concentration: 1.0 mg/ml

TREVIGEN[®]

8405 Helgerman Court, Gaithersburg, MD 20877 USA

Voice: 1-800-TREVIGEN (1-800-873-8443) • 301-216-2800

Fax: 301-216-2801 • e-mail: info@trevigen.com • www.trevigen.com

Streptavidin-AMCA

Catalog #: 4887-100-03

Storage: -20°C

TREVIGEN[®]

1-800-873-8443