

PRODUCT INSERT
**AFFINITY-ISOLATED ANTIBODIES
SPECIFIC TO GOAT IMMUNOGLOBULINS
UNCONJUGATED**

Product Code	Description	Volume	Antibody*
G50000	Swine Anti-Goat IgG (H+L)	2.0 ml	2.0 mg

Lot No.: See label **Expiration:** See label

Preparation:

Antibodies were affinity purified by column chromatography and adsorbed to remove cross reactivity to mouse and human immunoglobulins.

Quality Assurance:

Specificity and cross reactivities are tested in Caltag's in-house ELISA assay. Lot specific ELISA graphs are available upon request.

Preservatives:

This reagent contains *sodium azide*. Sodium azide yields hydrazoic acid under acid conditions, which is an extremely toxic compound. Azide compounds should be flushed with running water after being discarded. This procedure is recommended to

avoid deposits in plumbing where explosive conditions can develop.

Storage & Handling:

Store +2 to +8°C for short term (to 3 month) and -10°C or colder for long term. Aliquot and avoid repeated freezing and thawing. Mix thoroughly after thawing before use.

If the reagent is being diluted, it is recommended that only the quantity to be used within one week be diluted.

* Antibody value assigned is based on the extinction coefficient of IgG: Optical Density at 280 nm of 1.4 equals 1 mg antibody.

FOR RESEARCH USE ONLY. . . NOT FOR THERAPEUTIC OR IN VITRO DIAGNOSTIC USE

www.invitrogen.com

Invitrogen Corporation • 542 Flynn Rd • Camarillo • CA 93012 • Tel: 800.955.6288 • E-mail: techsupport@invitrogen.com

PI: L11058

(Rev 12/08) 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.