

**PRODUCT INSERT**

**AFFINITY-ISOLATED ANTIBODIES  
SPECIFIC TO HUMAN IMMUNOGLOBULINS  
ALKALINE PHOSPHATASE CONJUGATED**

Product Code	Description	Form	Volume
H10008	Goat Anti-Human IgG (Fc Sp.)	Alk. Phos.	1.0 ml
H10308	Goat Anti-Human IgG (H+L)	Alk. Phos.	1.0 ml
H14008	Goat Anti-Human IgA (α)	Alk. Phos.	1.0 ml
H15008	Goat Anti-Human IgM (μ)	Alk. Phos.	1.0 ml
H15708	Goat Anti-Human IgE (ε)	Alk. Phos.	1.0 ml
H16008	Goat Anti-Human Kappa (F+B)	Alk. Phos.	1.0 ml
H17008	Goat Anti-Human Ig's (Polyvalent)	Alk. Phos.	1.0 ml

**Lot No.:** See label

**Expiration:** See label

**Stabilizer:** A highly purified BSA has been added as a stabilizing agent

**Preparation:** Antibodies were affinity purified by column chromatography and adsorbed to remove unwanted specificities. The resulting antibody was conjugated to alkaline phosphatase (Alk. Phos.) according to the glutaraldehyde method.

**STORAGE & HANDLING**

Store reagents at 2-8°C. It is recommended that investigators dilute only the quantity to be used within one week.

**Buffer:** 50 mM TRIS, 1 mM MgCl<sub>2</sub>, 0.1% sodium azide, and 40% glycerol, v/v

**PRODUCT QUALITY CONTROL**

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

**Preservative:** 0.1% *sodium azide*. Sodium azide is an extremely toxic and dangerous compound particularly when combined with acids or metals. Solutions containing sodium azide should be disposed of properly.

**ANALYTE SPECIFIC REAGENT. ANALYTICAL AND PERFORMANCE CHARACTERISTICS ARE NOT ESTABLISHED.**

[www.invitrogen.com](http://www.invitrogen.com)

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

PI: L11059

(Rev 12/08) DCC-08-1818

© 2007 Invitrogen Corporation. All right reserved. These products may be covered by one or more Limited Use Label Licenses (see the Invitrogen Catalog or [www.invitrogen.com](http://www.invitrogen.com)). By use of these products you accept the terms and conditions of all applicable Limited Use Label Licenses.