

Goat anti-rabbit IgG (H&L) AgaroseCatalog number: 55015
Unit size: 10 mg**Product Details**

Storage Conditions	2-6°C and kept from light.
Expiration Date	12 months upon receiving
Concentration	Please see COA for detailed information
Formulation	1X PBS with 0.01% sodium azide

Unit Details

Unit	55015 (10 mg)
------	---------------

Antibody Properties

Species Reactivity	Rabbit
Class	Secondary
Clonality	Polyclonal
Host	Goat

Chemical Properties

Molecular Weight	~150000
------------------	---------

Biological Properties

Stabilizer	None
Preparation	Goat anti-rabbit IgG (H+L) is produced in goat with pooled total rabbit IgG, and affinity purified with rabbit IgG coupled beads.
Application	Antigen affinity chromatography

Applications

Agarose conjugated antibodies are designed for the rapid and efficient collection of antibodies and proteins, including recombinant fusion proteins, from a complex protein mixture (see protocol). Antibody affinity gels are useful for selection of immunoglobulins and immunoprecipitation of antigens from sera, antisera, ascitic fluid or culture fluid, bacterial and mammalian cell lysates and fusion protein preparations.