
Wheat Gliadin Polyclonal Antibody

(Catalog # A57785)

Background

Gliadin is a class of proteins present in wheat and several other cereals within the grass genus *Triticum*. Gliadins and glutenins are the two main components of the gluten fraction of the wheat seed. This gluten is found in products such as wheat flour. Gliadin is the water-soluble component of gluten, while glutenin is insoluble. Both of these proteins are necessary to allow bread to rise properly in baking. They are also examples of food-derived pathogenesis. People with gluten-sensitive enteropathy (the severe form of which is coeliac disease) are sensitive to alpha, beta, and gamma gliadins. Those with wheat-dependent (WD) exercise-induced anaphylaxis, WD urticaria and Baker's asthma are sensitive to omega-gliadins.

Description

Wheat Gliadin Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

Liquid. 0.03% Proclin 300, 50% Glycerol, 0.01M PBS, PH 7.4.

Specificity

Plant

Isotype

IgG

Purification

>95%, Protein G purified

Immunogen

Wheat Gliadin (Native Protein)

Storage

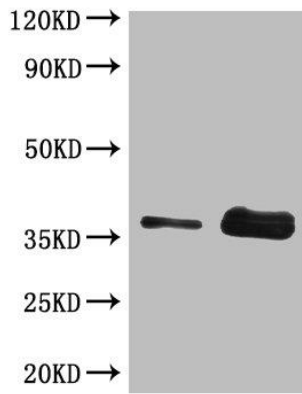
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

Wheat Gliadin

Application

ELISA, WB; Recommended dilution: WB:1:1000-1:5000



Western blot

All lanes: Wheat Gliadin Polyclonal Antibody at
2ug/ml+wheat flour

Lane 1: wheat flour at 2ug

Lane 2: wheat flour at 10ug

Secondary

Goat polyclonal to Rabbit IgG at 1/15000 dilution

Predicted band size: 35kDa

Observed band size: 35Da